

Agilent SLIMS 6.8

## Open Source Licenses

Agilent SLIMS contains software provided by third parties, including Open-Source software.  
This document contains licensing information about those third-party products used by Agilent SLIMS 6.8.

Title	Authors	Declared License
<b>AOP alliance</b>		public-domain
<b>Apache Commons BeanUtils</b>	Paul Jack, Stephen Colebourne, Berin Loritsch, Alex Crown, Marcus Zander, Paul Hamamnt, Rune Johannesen, Clebert Suconic, Norm Deane, Ralph Schaeer, Chris Audley, Rey François, Gregor Rayman, Jan Sorensen, Eric Pabst, Paulo Gaspar, Michael Smith, George Franciscus, Erik Meade, Tomas Viberg	Apache-2.0; Apache-1.1
<b>Apache Commons CLI</b>	Beluga Behr, Peter Donald, Brian Egge, Duncan Jones, Berin Loritsch, Peter Maddocks, Alexandru Mocanu, Andrew Shirley, Greg Thomas, Slawek Zachcial, Rubin Simons, James Strachan, Bob McWhirter, John Keyes, Rob Oxspring, Emmanuel Bourg, Thomas Neidhart, Rob Tompkins, Gary Gregory	Apache-2.0
<b>Apache Commons Codec</b>	Christopher O'Brien, Martin Redington, Jeffery Dever, Steve Zimmermann, Benjamin Walstrum, Oleg Kalnichevski, Dave Dribin, Alex Karasulu, Matthew Inger, Jochen Wiedmann, Sebastian Bazley, Matthew Pocock, Colm Rice, Adam Retter, Henri Yandell, Tim O'Brien, Scott Sanders, Rodney Waldhoff, Daniel Rall, Jon S. Stevens	Apache-2.0
<b>Apache Commons Codec</b>	Christopher O'Brien, Martin Redington, Jeffery Dever, Steve Zimmermann, Benjamin Walstrum, Oleg Kalnichevski, Dave Dribin, Alex Karasulu, Matthew Inger, Jochen Wiedmann, Sebastian Bazley, Matthew Pocock, Colm Rice, Adam Retter, Henri Yandell, Tim O'Brien, Scott Sanders, Rodney Waldhoff, Daniel Rall, Jon S. Stevens	Apache-2.0; Apache-1.1
<b>Apache Commons Compress</b>	Wolfgang Glas, Christian Kohlschütte, Bear Giles, Michael Kuss, Lasse Collin, John Kodis, BELUGA BEHR, Simon Spero, Michael Hausegger, Torsten Curdt, Stefan Bodewig, Sebastian Bazley, Christian Grobmeier, Julius Davies, Damjan Jovanovic, Emmanuel Bourg, Gary Gregory, Rob Tompkins, Peter Alfred Lee	Apache-2.0; Multi-license: Apache-2.0 OR bzip2-1.0.6; Multi-license: Apache-2.0 OR Info-ZIP
<b>Apache Commons CSV</b>	Henri Yandell, Martin van den Bemt, Yonik Seeley, Emmanuel Bourg, Gary Gregory, Benedikt Ritter, Rob Tompkins	Apache-2.0
<b>Apache Commons CSV</b>	Henri Yandell, Martin van den Bemt, Yonik Seeley, Emmanuel Bourg, Gary Gregory, Benedikt Ritter, Rob Tompkins	Apache-2.0; Apache-1.1
<b>Apache Commons DBCP</b>	Todd Carmichael, Wayne Woodfield, Dain Sundstrom, Philippe Mouawad, Glenn L. Nielsen, James House, James Ring, Peter Wicks, Morgan Delagrang, Geir Magnusson, Craig McClanahan, John McNally, Martin Poeschl, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Yoav Shapira, Jörg Schaible, Mark Thomas, Gary Gregory	Apache-2.0; Apache-1.1
<b>Apache Commons Exec</b>	Niklas Gustavsson, Benjamin Bentmann, Marco Ferrante, Jerome Lacoste, Milos Kleint, Pablo Hoertner, Niall Pemberton, Brett Porter, Trygve Laugstøl, Siegfried Goeschl, Sebastian Bazley, Gary Gregory	Apache-2.0; Apache-1.1
<b>Apache Commons FileUpload</b>	Aaron Freeman, Daniel Fabian, Jörg Heinicke, Stepan Koltsov, Michael Macaluso, Amichai Rothman, Alexander Sova, Paul Spurr, Thomas Vandahl, Henry Yandell, Jan Novotný, frank, maxxdev, Rafal Krzewski, Sean Legassick, Oleg Kalnichevski, David Sean Taylor, fangwentong, Martin Cooper, dlon Gillard	Apache-2.0; Apache-1.1
<b>Apache Commons IO</b>	Rahul Akolkar, Jason Anderson, Nathan Beyer, Emmanuel Bourg, Chris Eldredge, Magnus Grimsell, Jim Harrington, Thomas Ledoux, Andy Lehane, Marcelo Liberato, Alban Peignier, Adam Retter, Ian Springer, Dominik Stadler, Masato Tezuka, James Urie, Frank W. Zammetti, Scott Sanders, dlon Gillard, Nicola Ken Barozzi	Apache-2.0; Apache-1.1

<b>Apache Commons Lang</b>	C. Scott Ananian, Chris Audley, Stephane Bailliez, Michael Becke, Benjamin Bentmann, Ola Berg, Nathan Beyer, Stefan Bodewig, Janek Bogucki, Mike Bowler, Sean Brown, Alexander Day Chaffee, Al Chou, Greg Coladonato, Maarten Coene, Justin Couch, Michael Davey, Norm Deane, Morgan Delagrang, Ringo De Smet	Apache-2.0
<b>Apache Commons Logging</b>	Matthew P. Del Buono, Vince Eagen, Peter Lawrey, Berin Loritsch, Philippe Mouawad, Neeme Praks, Juozas Baliuka, Morgan Delagrang, Peter Donald, Robert Burrell Donkin, Simon Kitching, Dennis Lundberg, Costin Manolache, Craig McClanahan, Thomas Neidhart, Scott Sanders, Richard Sitze, Brian Stansberry, Rodney Waldhoff	Apache-2.0; Apache-1.1
<b>Apache Commons Validator</b>	Don Brown, Martin Cooper, David Graham, Ted Husted, Rob Leland, Craig McClanahan, James Mitchell, Niall Pemberton, James Turner, David Winterfeldt, Henri Yandell, Ben Speakmon, Nick Burch, Simone Tripodi, Benedikt Ritter, Gary Gregory	Apache-2.0
<b>Apache FontBox</b>		Apache-2.0; Multi-license: Apache-2.0 OR OFL-1.1; BSD-3-Clause
<b>Apache Groovy</b>	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0; BSD-3-Clause; MIT
<b>Apache Groovy</b>	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
<b>Apache Groovy</b>	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0; MIT; BSD-3-Clause
<b>Apache HttpClient</b>		Apache-2.0; Apache-1.1
<b>Apache HttpClient</b>		Apache-2.0; Apache-1.1
<b>Apache HttpClient Mime</b>		Apache-2.0; Apache-1.1
<b>Apache HttpCore</b>		Apache-2.0; Apache-1.1
<b>Apache Log4j</b>		Apache-2.0; Multi-license: Apache-2.0 OR TCL; Apache-1.1
<b>Apache Log4j API</b>		Apache-2.0; Apache-1.1
<b>Apache Log4j Core</b>		Apache-2.0; Multi-license: Apache-2.0 OR CDDL-1.0 OR MIT; Apache-1.1
<b>Apache Log4j SLF4J Binding</b>		Apache-2.0; Multi-license: Apache-2.0 OR MIT; Apache-1.1
<b>Apache Log4j Web</b>		Apache-2.0; Apache-1.1
<b>Apache Lucene (module: core)</b>		Multi-license: Apache-2.0 OR BSD-2-Clause OR BSD-3-Clause OR ICU OR MIT; Multi-license: Apache-2.0 OR MIT; Apache-2.0; BSD-3-Clause; MIT; ICU
<b>Apache Lucene (module: queryparser)</b>		Apache-2.0; Multi-license: Apache-2.0 OR MIT; Multi-license: Apache-2.0 OR BSD-2-Clause OR BSD-3-Clause OR ICU OR MIT
<b>Apache PDFBox</b>		Unicode; Zlib; BSD-3-Clause; APAFML; OFL-1.1; Apache-2.0

Apache PDFBox tools		Apache-2.0; Apache-1.1
Apache POI		Apache-2.0
Apache POI - API based on OPC and OOXML schemas		Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR EPL-1.0 OR MIT
Apache POI - Common		Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR EPL-1.0 OR MIT; MIT
Apache Tika parser modules		Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR CDDL-1.0 OR MIT OR MPL-1.1 OR NetCDF OR SAX-PD; Multi-license: Apache-2.0 OR ICU; Multi-license: Apache-2.0 OR MIT; Multi-license: Apache-2.0 OR IPA; Apache-1.1
Apache Velocity	Will Glass-Husain, Geir Magnusson Jr., Daniel Rall, Henning P. Schmiedehausen, Nathan Bubna	Apache-2.0; Apache-1.1
Apache XBean :: Spring		Apache-2.0
Apache XML Security for Java	Sean Mullan, Berin Lautenbach, Axl Mattheus, Erwin van der Koog, Raul Benito Garcia, Vishal Mahajan, Milan Tomic, Colm O hEigeartaigh	Apache-2.0; Multi-license: Apache-2.0 OR MIT; W3C-19980720; Apache-1.1; W3C
ARQ		BSD-2-Clause; BSD-3-Clause; JSON
ASM based accessors helper used by json-smart	Uriel Chemouni, Zhang Jian He	Apache-2.0
AutoValue Annotations		Apache-2.0
barbecue	Sean C. Sullivan, Werner Keil	BSD-3-Clause
Barcode4J	Jeremias Märki, Web Maestro Clay Leeds, Nicola Ken Barozzi, Dietmar Bürkle	Apache-2.0
batik-awt-util		Apache-2.0
batik-bridge		Apache-2.0
batik-dom		Apache-2.0; W3C
batik-svggen		Apache-2.0
batik-util		Apache-2.0
batik-xml		Apache-2.0
Bean Validation API	Emmanuel Bernard, Hardy Ferentschik, Gunnar Morling, Guillaume Smet	Apache-2.0
Bouncy Castle Provider		BouncyCastle; MIT
Bouncy Castle S/MIME API		
Bouncy Castle TSP/MIME API		BouncyCastle
Camel :: Attachments		Apache-2.0
Camel :: Core		Apache-2.0
Camel :: HazelCast		Apache-2.0
Camel :: JMS		Apache-2.0
Camel :: Mail		Apache-2.0
Camel :: Management		Apache-2.0
Camel :: Netty		Apache-2.0
Camel :: Quartz		Apache-2.0
Camel :: Spring		Apache-2.0
Camel :: Spring Integration		Apache-2.0
Camel :: Spring Main		Apache-2.0
Camel :: Support		Apache-2.0

Camel :: Test (d eprecated)		Apache-2.0
Camel :: Test :: Spring (d eprecated)		Apache-2.0
cglib		Multi-license: Apache-2.0 OR unknown; Apache-2.0; BSD-3-Clause
CGLIB	Juozas Baliuka, Chris Nokleberg, Matas Ramoska, Sam Berlin, Wes Biggs	Apache-2.0; BSD-3-Clause
Commons DBCP	Todd Carmichael, Wayne Woodfield, Dain Sundstrom, Philippe Mouawad, Morgan Delagrang, Geir Magnusson, Craig McClanahan, John McNally, Martin Poeschl, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Yoav Shapira, Phil Steitz, Jörg Schaible, Mark Thomas	Apache-2.0
Commons Discovery	Richard Sitze, Craig R. McClanahan, Costin Manolache, James Strachan, Matthew Hawthorne, Davanum Srinivas, Rory Winston, Robert Burrell Donkin, Simone Tripodi	Apache-2.0
Commons JXPath	Dmitri Plotnikov, Craig McClanahan, Matt Benson	Apache-2.0
Commons Lang	C. Scott Ananian, Chris Audley, Stephane Bailliez, Michael Becke, Benjamin Bentmann, Ola Berg, Nathan Beyer, Stefan Bodewig, Janek Bogucki, Mike Bowler, Sean Brown, Alexander Day Chaffee, Al Chou, Greg Coladonato, Maarten Coene, Justin Couch, Michael Davey, Norm Deane, Ringo De Smet, Russel Dittmar	Apache-2.0
Commons Math	C. Scott Ananian, Mark Anderson, Rémi Amtzen, Michael Bjorkegren, John Bollinger, Cyril Briquet, Paul Cowan, Benjamin Croizet, Larry Diamond, Rodrigo di Lorenzo Lopes, Hasan Diwan, Ted Dunning, John Gant, Ken Geis, Bernhard Grünewaldt, Elliotte Rusty Harold, Matthias Hummel, Ismael Juma, Eugene Kirpichov, Piotr Kochanski	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause; Multi-license: Apache-2.0 OR BSD-2-Clause
Date Calculator Common		Apache-2.0
Date Calculator for Joda		Apache-2.0
dom4j	Curt Arnold, David Lucas, David White, Jakob Jenkov, James Strachan, Laramie Crocker, Maarten Coene, Michael Skells, Nick Sanderson, Steen Lehmann, Tobias Rademacher, Bob McWhirter, Todd Wolff, OuYang Chen, Yuxin Ruan	dom4j; proprietary-license
domino-jackson		Apache-2.0
domino-jackson-super		Apache-2.0
Elemental2 Core		Apache-2.0
Elemental2 Dom		Apache-2.0
Elemental2 Promise		Apache-2.0
EventBinder		Apache-2.0
FindBugs-jsr305		Apache-2.0; CC-BY-2.5
fr.opensagres.xdocreport.converter.docx.xwpf		MIT
GIN (GWT INjection)	Brian Stoler, Peter Schmitt, Jesse Wilson, Robbie Vanbrabant, Fazal Asim, Daniel Burrows, Scott Blum, Adam Winer, Philippe Beaudoin	Apache-2.0
Google Guice - Core Library		Multi-license: Apache-2.0 OR CC0-1.0; Apache-1.1; Apache-2.0
Google Guice - Extensions - AssistedInject		Apache-1.1; Apache-2.0
Guava: Google Core Libraries for Java		CC0-1.0; Apache-2.0
Guava GWT compatible libs		Apache-2.0; Multi-license: Apache-2.0 OR CC0-1.0; CC0-1.0
GWT CLDR		Apache-2.0

GWT DateTimeFormat		Apache-2.0
GWT i18n		Apache-2.0
GWT Servlet Library 1.5		Apache-2.0; CC0-1.0
gwt-user		Apache-2.0; public-domain
Hamcrest All		BSD-3-Clause-Clear
HAPI - 01 - Base Module		MPL-1.1; Zlib
HAPI - Java HL7 API - Generated Structures v2.7		Multi-license: GPL-1.0-or-later OR MPL-1.1; Multi-license: GPL-1.0-or-later OR MPL-1.1 OR OML; OML
HAPI - The Open Source Java HL7 Parser and API	Axel Biernat, Rodolfo Chavez, Ryan W. Gross, Nicolas Lefebvre, Jens Kristian Villadsen, Bryan Tripp, James Agnew, Christian Ohr, Niranjana Sharma, Alexei Guevara, Charles D. Fisher, Abraham Tio, Leslie Mann, Neal Acharya, Nico Vannieuwenhuyze	Multi-license: GPL-3.0-only OR MPL-1.1 OR MPL-2.0; MPL-1.1
hazelcast		Apache-2.0; MIT; AML; CC0-1.0
hazelcast-spring		Apache-2.0; BSD-3-Clause
Hibernate Commons Annotations		LGPL-2.1-or-later
hibernate-core-relocation		Multi-license: LGPL-2.1-only OR LGPL-2.1-or-later; LGPL-2.1-or-later; Multi-license: BSD-3-Clause OR EPL-2.0; CC0-1.0
hibernate-core-relocation		LGPL-3.0-only
Hibernate Validator Engine - Relocation Artifact		Apache-2.0
Hibernate Validator Engine - Relocation Artifact		Apache-2.0
HttpClient	Armando Anton, Sebastian Bazley, Ola Berg, Sam Berlin, Mike Bowler, Samit Jain, Eric Johnson, Christian Kohlschuetter, Ryan Lubke, Sam Maloney, Rob Di Marco, Juergen Pill, Mohammad Rezaei, Roland Weber, Laura Werner, Mikael Wilstrom, Michael Becke, Jeff Dever, dlon Gillard, Ortwin Glueck	Apache-2.0; Multi-license: Apache-2.0 OR MIT
ICU4J	Mark Davis, John Emmons, Doug Felt, Deborah Goldsmith, Steven Loomis, Markus Scherer, Peter Edberg, Yoshito Umaoka	Multi-license: BSD-2-Clause OR BSD-3-Clause OR public-domain OR X11; Unicode-DFS-2016; Multi-license: Unicode-DFS-2016 OR X11; Multi-license: IPA OR Unicode-DFS-2016 OR X11
Jackson-annotations		Apache-2.0
Jackson-core		Apache-2.0
jackson-databind		Apache-2.0
Jackson-dataformat-CSV		Apache-2.0
Jackson-dataformat-YAML		Apache-2.0
Jackson-dataformat-YAML		Apache-2.0
Jackson datatype: Guava		Apache-2.0
Jackson datatype: jdk8		Apache-2.0
Jackson datatype: JSR310		Apache-2.0
JasperReports Font Extension		LGPL-3.0-only
JasperReports Functions		GPL-3.0-or-later
JasperReports Library	Teodor Danciu, Lucian Chirita	LGPL-3.0-only; GPL-3.0-or-later; LGPL-3.0-or-later
Java Annotation Indexer		Apache-2.0

<b>JavaBeans(TM) Activation Framework</b>		CDDL-1.0; Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception
<b>java.jwt</b>	Auth0, Hernan Zalazar, Luciano Balmaceda	MIT
<b>JavaMail API</b>		Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.0 OR CDDL-1.1 OR GPL-2.0-with-classpath-exception; Apache-2.0
<b>Java Native Access</b>	Timothy Wall, Matthias Bläsing	Multi-license: Apache-2.0 OR LGPL-2.1-or-later
<b>Java Persistence API, Version 2.1</b>	Emmanuel Bernard, Hardy Ferentschik, Steve Ebersole	EPL-1.0; Multi-license: BSD-3-Clause OR EPL-1.0
<b>Java Servlet API</b>	Ed Burns, Shing Wai Chan	Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Apache-2.0; Multi-license: Apache-2.0 OR CDDL-1.1 OR GPL-2.0-with-classpath-exception; Multi-license: Apache-2.0 OR CDDL-1.1 OR GPL-2.0-with-classpath-exception OR IETF
<b>Java Servlet API</b>	Ed Burns, Shing Wai Chan	Multi-license: CDDL-1.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Multi-license: Apache-2.0 OR CDDL-1.1 OR GPL-2.0-with-classpath-exception
<b>Javassist</b>	Shigeru Chiba, Andrew Dinn, Kabir Khan, Scott Marlow	
<b>javax.annotation-api</b>		Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Apache-2.0
<b>jaxb-api</b>	Martin Grebac, Iaroslav Savitskiy	Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Apache-2.0
<b>jaxb-api</b>		
<b>jaxen</b>	Ryan Gustafson, David Peterson, Mark Wilson, Jacob Kjome, Michael Brennan, Jason Hunter, Brett McLaughlin, Bradley S. Huffman, K. Ari Krupnikov, Paul R. Brown, Guoliang Cao, Jérôme Nègre, Eddie McGreal, Steen Lehmann, Ben McCann, Don Corley, Bob McWhirter, James Strachan, David Megginson, Erwin Bolwidt	BSD-3-Clause; Apache-2.0
<b>jCIFS</b>		LGPL-2.1-only; LGPL-2.1-or-later; BSD-2-Clause; LGPL-2.0-or-later
<b>jCSV</b>		Apache-2.0
<b>JDOM</b>	Jason Hunter, Rolf Lear	Apache-2.0
<b>Jena</b>		BSD-2-Clause; BSD-3-Clause; AML
<b>JGraphT - Core</b>		Multi-license: EPL-1.0 OR LGPL-2.1-only; EPL-1.0
<b>Joda-Time</b>	Adrian Acala, Afif Ahmed, AlexMe951, Guy Allard, Brad Arndt, Mikey Austin, Oren Benjamin, Fredrik Borgh, Dave Brosius, Dan Cavallaro, Luc Claes, Emiliano Claria, Dan Cojocar, Evgeniy Devyatkykh, dspitfire, Christopher Elkins, emopers, Michael Ernst, Jeroen van Erp, Gwyn Evans	Apache-2.0; Multi-license: Apache-2.0 OR MIT
<b>JsInterop Annotations</b>		Apache-2.0

Js Interop Base		Apache-2.0
JSON.simple		Apache-2.0
JSON Small and Fast Parser	Uriel Chemouni, Eitan Raviv, ZhangJian He	Apache-2.0
jsoup Java HTML Parser		MIT
JTA 1.1		Apache-2.0
JUnit	David Saff, Kevin Cooney, Stefan Birkner, Marc Philipp	EPL-1.0
Kotlin Stdlib Jdk8		Apache-2.0
Kubernetes Discovery Plugin for Hazelcast		Apache-2.0
Liquibase		Apache-2.0; Multi-license: Apache-2.0 OR BSD-2-Clause OR CDDL-1.0 OR CDDL-1.1 OR CPL-1.0 OR ISC OR LGPL-3.0-only OR MIT OR Plexus; LGPL-3.0-only; CDDL-1.0; MIT; Plexus; CPL-1.0; ISC; BSD-2-Clause
liquibase-slf4j		MIT
Lucene Common Analyzers		Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause; Multi-license: Apache-2.0 OR BSD-2-Clause OR BSD-3-Clause OR ICU OR MIT; Multi-license: Apache-2.0 OR MIT; Multi-license: Apache-2.0 OR IPA; BSD-3-Clause
mysql-java-client	Vladislav Vaintroub, Massimo Siani, Georg Richter, Diego Dupin	Multi-license: BSD-3-Clause OR LGPL-2.1-only OR LGPL-2.1-or-later; Multi-license: BSD-3-Clause OR LGPL-2.1-or-later
micrometer-core		Apache-2.0; Multi-license: Apache-2.0 OR MIT
micrometer-registry-new-relic		Apache-2.0
micrometer-registry-prometheus		Apache-2.0
Microsoft JDBC Driver for SQL Server		MIT; Apache-2.0
mockito-core	Szczepan Faber, Brice Dutheil, Rafael Winterhalter, Tim van der Lippe	MIT; Apache-2.0; BSD-3-Clause
mockito-core	Szczepan Faber, Brice Dutheil, Rafael Winterhalter, Tim van der Lippe	MIT; Apache-2.0
Netty/All-in-One		Apache-2.0
OAuth2 for Spring Security		Apache-2.0
Objenesis		Apache-2.0
Old JAXB Core		Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Apache-2.0; Multi-license: CDDL-1.0 OR CDDL-1.1 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception OR SAX-PD
Old JAXB Runtime		Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception OR public-domain
org.apache.poi.xwpf.converter.pdf		MIT
OWASP Java HTML Sanitizer		BSD-3-Clause



quartz		Apache-2.0
REngine Java interface to R		Multi-license: GPL-2.0-only OR LGPL-2.1-only
Rserve client back-end for REngine		Multi-license: GPL-2.0-only OR LGPL-2.1-only
Spring Beans		Apache-2.0
spring-boot-starter-thymeleaf		
spring-boot-starter-web		
Spring Context		Apache-2.0
Spring Core		Apache-2.0; BSD-3-Clause
springfox-core	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-swagger2	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-swagger-ui	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
Spring HATEOAS	Oliver Drotbohm, Greg Turnquist	Apache-2.0
Spring LDAP	Eric Dalquist, Marius Scurtescu, Tim Terry, Keith Barlow, Paul Harvey, Marvin S. Addison, Mattias Hellborg, Arthursson, Ulrik Sandberg	Apache-2.0
Spring Messaging		Apache-2.0
Spring Object/Relational Mapping		Apache-2.0
spring-security-config	Rob Winch, Joe Grandja	Apache-2.0
spring-security-core	Rob Winch, Joe Grandja	Apache-2.0
spring-security-ldap	Rob Winch, Joe Grandja	Apache-2.0
spring-security-oauth2-client	Rob Winch, Joe Grandja	Apache-2.0
spring-security-oauth2-jose	Rob Winch, Joe Grandja	Apache-2.0
spring-security-saml2-service-provider	Rob Winch, Joe Grandja	Apache-2.0
spring-security-web	Rob Winch, Joe Grandja	Apache-2.0
Spring TestContext Framework		Apache-2.0
Spring Transaction		Apache-2.0
Spring Web		Apache-2.0
Spring Web MVC		Apache-2.0
swagger-annotations		Apache-2.0
swagger-models		
thymeleaf	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0
thymeleaf-spring5	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0
vaadin-server	Artur Signell, Marko Gronroos	Apache-2.0
vaadin-spring		Apache-2.0
vaadin-themes	Artur Signell, Marko Gronroos	Apache-2.0
VT Password Libraries	Daniel Fisher, Marvin Addison	Apache-2.0
Xalan Java		W3C-19980720; Multi-license: Apache-2.0 OR TCL; Apache-2.0
Xerces2-j		W3C-19980720; Apache-2.0; Multi-license: Apache-2.0 OR IPA
ZXing Core		Apache-2.0
Deep Dependencies		
Title	Authors	Declared License
Adobe XMPCore		BSD-2-Clause
AntLR		BSD-2-Clause

<b>Apache Ant + ANTLR</b>		Apache-2.0; Multi-license: Apache-2.0 OR SAX-PD OR W3C; Apache-1.1
<b>Apache Ant Core</b>		Multi-license: Apache-2.0 OR SAX-PD OR W3C; Apache-2.0; Info-ZIP
<b>Apache Ant + JUnit</b>		Apache-2.0; Multi-license: Apache-2.0 OR SAX-PD OR W3C; Apache-1.1
<b>Apache Ant Launcher</b>		Apache-2.0; Multi-license: Apache-2.0 OR SAX-PD OR W3C
<b>Apache Commons BeanUtils</b>	Paul Jack, Stephen Colebourne, Berin Loritsch, Alex Crown, Marcus Zander, Paul Hamamnt, Rune Johannesen, Clebert Suconic, Norm Deane, Ralph Schaer, Chris Audley, Rey François, Gregor Rayman, Jan Sorensen, Eric Pabst, Paulo Gaspar, Michael Smith, George Franciscus, Erik Meade, Tomas Viberg	Apache-2.0; Apache-1.1
<b>Apache Commons Codec</b>	Christopher O'Brien, Martin Redington, Jeffery Dever, Steve Zimmermann, Benjamin Walstrum, Oleg Kalnichevski, Dave Dribin, Alex Karasulu, Matthew Inger, Jochen Wiedmann, Sebastian Bazley, Matthew Pocock, Colm Rice, Adam Retter, Henri Yandell, Tim O'Brien, Scott Sanders, Rodney Waldhoff, Daniel Rall, Jon S. Stevens	Apache-2.0; Apache-1.1
<b>Apache Commons Collections</b>	Rafael U. C. Afonso, Max Rydahl Andersen, Avalon, Federico Barbieri, Jeffrey Barnes, Nicola Ken Barozzi, Arron Bates, Sebastian Bazley, Benjamin Bentmann, Ola Berg, Sam Berlin, Christopher Berry, Nathan Beyer, Rune Peter Bjørnstad, Janek Bogucki, Maarten Brak, Dave Bryson, Chuck Burdick, Julien Buret, Josh Cain	Apache-2.0
<b>Apache Commons Collections</b>	Rafael U. C. Afonso, Max Rydahl Andersen, Avalon, Federico Barbieri, Jeffrey Barnes, Nicola Ken Barozzi, Arron Bates, Sebastian Bazley, Benjamin Bentmann, Ola Berg, Sam Berlin, Christopher Berry, Nathan Beyer, Rune Peter Bjørnstad, Janek Bogucki, Maarten Brak, Dave Bryson, Chuck Burdick, Julien Buret, Josh Cain	Apache-2.0; Apache-1.1
<b>Apache Commons Compress</b>	Wolfgang Glas, Christian Kohlschütte, Bear Giles, Michael Kuss, Lasse Collin, John Kodis, BELUGA BEHR, Simon Spero, Michael Hausegger, Torsten Curdt, Stefan Bodewig, Sebastian Bazley, Christian Grobmeier, Julius Davies, Damjan Jovanovic, Emmanuel Bourg, Gary Gregory, Rob Tompkins, Peter Alfred Lee	Apache-2.0; Multi-license: Apache-2.0 OR Info-ZIP; Apache-1.1
<b>Apache Commons IO</b>	Rahul Akolkar, Jason Anderson, Nathan Beyer, Emmanuel Bourg, Chris Eldredge, Magnus Grimsell, Jim Harrington, Thomas Ledoux, Andy Lehane, Marcelo Liberato, Alban Peignier, Adam Retter, Ian Springer, Dominik Stadler, Masato Tezuka, James Urie, Frank W. Zammetti, Scott Sanders, Dion Gillard, Nicola Ken Barozzi	Apache-2.0; Apache-1.1
<b>Apache Commons Lang</b>	C. Scott Ananian, Chris Audley, Stephane Bailliez, Michael Becke, Benjamin Bentmann, Ola Berg, Nathan Beyer, Stefan Bodewig, Janek Bogucki, Mike Bowler, Sean Brown, Alexander Day Chaffee, Al Chou, Greg Coladonato, Maarten Coene, Justin Couch, Michael Davey, Norm Deane, Morgan Delagrang, Ringo De Smet	Apache-2.0; Apache-1.1
<b>Apache Commons Math</b>	Eldar Agalarov, Tim Allison, C. Scott Ananian, Mark Anderson, Peter Andrews, Rémi Arntzen, Matt Adereth, Jared Becksfort, Michael Bjorkegren, Brian Bloniarz, John Bollinger, Cyril Briquet, Dave Brosius, Dan Checkoway, Anders Conbere, Charles Cooper, Paul Cowan, Benjamin Croizet, Larry Diamond, Aleksei Dievskii	Multi-license: Apache-2.0 OR BSD-3-Clause; Apache-2.0; BSD-3-Clause; BSD-2-Clause; Apache-1.1
<b>Apache Commons Pool</b>	Morgan Delagrang, Geir Magnusson, Craig McClanahan, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Robert Burrell Donkin, Sandy McArthur, Simone Tripodi, Gary Gregory, Matt Sicker	Apache-2.0; Apache-1.1; Multi-license: Apache-2.0 OR cc-pd
<b>Apache Geronimo JMS Spec 2.0</b>		Apache-2.0; Apache-1.1



Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0

Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache Groovy	Joern Eyrich, Robert Kuzelj, Rod Cope, Yuri Schimke, James Birchfield, Robert Fuller, Sergey Udovenko, Hallvard Traetteberg, Peter Reilly, Brian McCallister, Richard Monson-Haefel, Brian Larson, Artur Biesiadowski, Ivan Z. Ganza, Larry Jacobson, Jake Gage, Arjun Nayyar, Masato Nagai, Mark Chu-Carroll, Mark Turansky	Apache-2.0
Apache HttpClient		Apache-2.0; Apache-1.1
Apache HttpCore		Apache-2.0; Apache-1.1
Apache JempBox		Apache-2.0
Apache Lucene (module: queries)		Apache-2.0; Multi-license: Apache-2.0 OR BSD-2-Clause OR BSD-3-Clause OR ICU OR MIT; Multi-license: Apache-2.0 OR MIT
Apache Lucene (module: sandbox)		Apache-2.0; Multi-license: Apache-2.0 OR BSD-2-Clause OR BSD-3-Clause OR ICU OR MIT; Multi-license: Apache-2.0 OR MIT
Apache PDFBox Debugger		Apache-2.0; Apache-1.1
Apache POI		Multi-license: Apache-2.0 OR BSD-3-Clause OR EPL-1.0 OR MIT; Apache-2.0
Apache Preflight		Apache-2.0; Apache-1.1
Apache Tika core		Apache-2.0; Apache-1.1
Apache XML Graphics Commons		Apache-2.0
Apache XmpBox		Apache-2.0; Apache-1.1
asm	Eric Bruneton, Eugene Kuleshov, Remi Forax	BSD-3-Clause; Apache-2.0
AssertJ fluent assertions		Apache-2.0
attoparser		Apache-2.0
Avalon Framework Implementation		Apache-2.0
batik-anim		Apache-2.0
batik-constants		Apache-2.0
batik-css		Apache-2.0
batik-ext		Apache-2.0; W3C
batik-gvt		Apache-2.0
batik-i18n		Apache-2.0
batik-parser		Apache-2.0
batik-script		Apache-2.0; Multi-license: Apache-2.0 OR TCL
batik-shared-resources		Apache-2.0
batik-svg-dom		Apache-2.0; MIT; W3C

Bouncy Castle ASN.1 Extension and Utility APIs		
Bouncy Castle PKIX, CMS, EAC, TSP, PKCS, OCSP, CMP, and CRMF APIs		
Byte Buddy (without dependencies)		BSD-3-Clause; Apache-2.0
c3p0		Multi-license: LGPL-2.1-only OR EPL-1.0; LGPL-2.1-only
Camel::API		Apache-2.0
Camel::Base		Apache-2.0
Camel::Base Engine		Apache-2.0
Camel::Bean		Apache-2.0
Camel::Browse		Apache-2.0
Camel::Cluster		Apache-2.0
Camel::Controlbus		Apache-2.0
Camel::Core Engine		Apache-2.0
Camel::Core Languages		Apache-2.0
Camel::Core Model		Apache-2.0
Camel::Core Processor		Apache-2.0
Camel::Core Reifier		Apache-2.0
Camel::Core XML		Apache-2.0
Camel::Dataformat		Apache-2.0
Camel::DataSet		Apache-2.0
Camel::Direct		Apache-2.0
Camel::Direct VM		Apache-2.0
Camel::File		Apache-2.0
Camel::Health		Apache-2.0
Camel::Language		Apache-2.0
Camel::Log		Apache-2.0
Camel::Main		Apache-2.0
Camel::Management API		Apache-2.0
Camel::Mock		Apache-2.0
Camel::Ref		Apache-2.0
Camel::REST		Apache-2.0
Camel::Saga		Apache-2.0
Camel::Scheduler		Apache-2.0
Camel::Seda		Apache-2.0
Camel::SpringXML		Apache-2.0
Camel::Stub		Apache-2.0
Camel::Timer		Apache-2.0
Camel::Tooling::Model		Apache-2.0
Camel::Util		Apache-2.0
Camel::Util::JSon		Apache-2.0
Camel::Validator		Apache-2.0
Camel::VM		Apache-2.0

Camel::XML IO Util		Apache-2.0
Camel::XML JAXB		Apache-2.0
Camel::XML JAXP		Apache-2.0
Camel::XPath		Apache-2.0
Camel::XSLT		Apache-2.0
Castor CORE - Core code/functionality		Apache-2.0; dom4j
Castor XML - core		dom4j; Apache-2.0
Checker Qual	Michael Ernst, Suzanne Millstein	MIT
Checker Qual	Michael Ernst, Suzanne Millstein	MIT
ClassMate	Tatu Saloranta, Brian Langel	Apache-2.0
ClassMate	Tatu Saloranta, Brian Langel	Apache-2.0
commons-collections	Max Rydahl Andersen, Federico Barbieri, Arron Bates, Nicola Ken Barozzi, Ola Berg, Christopher Berry, Janek Bogucki, Chuck Burdick, Dave Bryson, Julien Buret, Jonathan Carlson, Ram Chidambaram, Peter Donald, Steve Downey, Rich Dougherty, Stefano Fornari, Andrew Freeman, Gerhard Froehlich, Paul Jack, Eric Johnson	Apache-2.0
Commons Digester	Bradley M. Handy, Christopher Lenz, Ted Husted, David H. Martin, Henri Chen, Janek Bogucki, Mark Huisman, Paul Jack, Anton Maslovsky, Matt Cleveland, Gabriele Carcassi, Wendy Smoak, Kevin Ross, Craig McClanahan, Robert Burrell Donkin, Scott Sanders, James Strachan, Jason van Zyl, Tim O'Brien, Jean-Francois Arcand	Apache-2.0
Commons Digester	Bradley M. Handy, Christopher Lenz, Ted Husted, David H. Martin, Henri Chen, Janek Bogucki, Mark Huisman, Paul Jack, Anton Maslovsky, Matt Cleveland, Gabriele Carcassi, Wendy Smoak, Kevin Ross, Craig McClanahan, Robert Burrell Donkin, Scott Sanders, James Strachan, Jason van Zyl, Tim O'Brien, Jean-Francois Arcand	Apache-2.0
Commons Pool	Morgan Delagrang, Geir Magnusson, Craig McClanahan, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Robert Burrell Donkin, Sandy McArthur, Gary Gregory	Apache-2.0
Commons Pool	Morgan Delagrang, Geir Magnusson, Craig McClanahan, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Robert Burrell Donkin, Sandy McArthur, Gary Gregory	Apache-2.0
Cryptacular Library	Daniel Fisher, Marvin S. Addison	Multi-license: Apache-2.0 OR LGPL-3.0-only
curvesapi	stormdollar, Dustin Spicuzza	BSD-2-Clause; BSD-3-Clause; Apache-2.0
dec		MIT
dom4j		dom4j
Eclipse Compiler for Java(TM)		Multi-license: EPL-1.0 OR EPL-2.0; EPL-2.0; EPL-1.0
Elemental2 Promise		Apache-2.0
error-prone annotations		Apache-2.0
Extended StAX API	Roman Grigoriadi, Zheng Jun Li	Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Apache-2.0
fastinfoset		Apache-2.0; Multi-license: Apache-1.1 OR Apache-2.0; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception
Flute		W3C; W3C-19980720
fr.opensagres.poi.xwpf.converter.core		MIT

fr.opensagres.poi.xwpf.converter.pdf		MIT
fr.opensagres.poi.xwpf.converter.html		MIT
fr.opensagres.xdocreport.converter		MIT
fr.opensagres.xdocreport.core		MIT; Multi-license: Apache-2.0 OR MIT
fr.opensagres.xdocreport.itext.extension		MIT
GenTyRef		Apache-2.0
Guava: Google Core Libraries for Java		Apache-2.0; CC0-1.0
Guava InternalFutureFailureAccess and InternalFutures		Apache-2.0
Guava ListenableFuture only		Apache-2.0
gwt-elemental		Apache-2.0
Hamcrest Core		BSD-3-Clause-Clear
HdrHistogram		Multi-license: BSD-2-Clause OR CC0-1.0; CC0-1.0
Hibernate Commons Annotations		
HikariCP		Apache-2.0
ICU4J	Mark Davis, John Emmons, Doug Felt, Deborah Goldsmith, Steven Loomis, Markus Scherer, Peter Edberg, Yoshito Umaoka	ICU; MIT; GPL-3.0-only
ISO Parser		Apache-2.0; EPL-1.0
istack common utility code runtime		Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Apache-2.0; Multi-license: CDDL-1.0 OR CDDL-1.1 OR GPL-2.0-with-classpath-exception; CDDL-1.1
J2ObjC Annotations		Apache-2.0
Jackson-annotations		Apache-2.0
Jackson BOM		Apache-2.0
Jackson-core		Apache-2.0
jackson-databind		Apache-2.0
Jakarta Activation		BSD-3-Clause; Multi-license: EPL-2.0 OR GPL-2.0-with-classpath-exception
Jakarta Activation API		BSD-3-Clause; Multi-license: EPL-2.0 OR GPL-2.0-with-classpath-exception
Jakarta Mail API		BSD-3-Clause; Multi-license: EPL-2.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception
jakarta-regexp		Apache-2.0
Jakarta XML Binding API		BSD-3-Clause; Multi-license: Apache-2.0 OR BSD-3-Clause OR MIT; Multi-license: EPL-2.0 OR GPL-2.0-with-classpath-exception
Java Advanced Imaging Image I/O Tools API core (standalone)	John Hewson, Peter Hull, Mark Carroll, Robin Stevens, Yannick De Turck, Butch Howard	BSD-2-Clause; BSD-3-Clause-No-Nuclear-Warranty
Java Annotation Indexer		Apache-2.0
JavaBeans Activation Framework		Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-



		2.0-with-classpath-exception; Apache-2.0
JavaBeans Activation Framework API jar		Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Apache-2.0
JavaBeans(TM) Activation Framework		CDDL-1.0
javaparser-core		Multi-license: Apache-2.0 OR LGPL-3.0-only
JavaServlet(TM) Specification		Apache-2.0
Javassist	Shigeru Chiba, Andrew Dinn, Kabir Khan, Scott Marlow	Multi-license: Apache-2.0 OR LGPL-2.1-or-later OR MPL-1.1
java-support		Apache-2.0
javax.inject		Apache-2.0
javax.persistence-api		Multi-license: BSD-3-Clause OR EPL-1.0; Apache-2.0
JAXB Runtime		CDDL-1.0; Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception OR SAX-PD; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; CDDL-1.1; Apache-2.0
JBoss Logging3		Apache-2.0
jboss-transaction-api_1.2_spec		
JCIP Annotations under Apache License		Apache-2.0
JCL 1.2 implemented over SLF4J		Apache-2.0; MIT
JCL 1.2 implemented over SLF4J		Apache-2.0
jcommander		Apache-2.0
JCommon		Multi-license: LGPL-2.1-or-later OR LGPL-3.0-only; LGPL-2.1-or-later; Apache-2.0; LGPL-2.1-only; LGPL-2.0-or-later; GPL-2.0-only
Jena IRI		BSD-2-Clause; BSD-3-Clause
JetBrains Java Annotations		Apache-2.0
JFreeChart		LGPL-3.0-only; LGPL-2.1-or-later; BSD-3-Clause
JLine	Scott Gray, Hiram Chirino, Jason Dillon, Guillaume Nodet, Marc Prud'hommeaux	BSD-3-Clause; Apache-2.0
JMatIO		BSD-2-Clause
Joda-Time	Adrian Acala, Afif Ahmed, AlexMe951, Guy Allard, Brad Arndt, Mikey Austin, Oren Benjamin, Fredrik Borgh, Dave Brosius, Dan Cavallaro, Luc Claes, Emiliano Claria, Dan Cojocar, Evgeniy Devyatykh, dspitfire, Christopher Elkins, emopers, Michael Ernst, Jeroen van Erp, Gwyn Evans	Apache-2.0
jquery		MIT
JsInterop Annotations		Apache-2.0
JsInterop Base		Apache-2.0
JTidy		W3C-19980720; MIT
JUL to SLF4J bridge		MIT
JUL to SLF4J bridge		MIT
JUnit	David Saff, Kevin Cooney, Stefan Birkner, Marc Philipp	EPL-1.0

<b>JUnit 5 (Bill of Materials)</b>	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0
<b>JUnit Jupiter API</b>	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-2.0
<b>JUnit Jupiter Engine</b>	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-2.0
<b>JUnit Platform Commons</b>	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-2.0
<b>JUnit Platform Engine API</b>	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-2.0
<b>JUnit Platform Launcher</b>	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0
<b>Kotlin Stdlib</b>		Apache-2.0
<b>Kotlin Stdlib Common</b>		Apache-2.0
<b>Kotlin Stdlib Jdk7</b>		Apache-2.0
<b>LatencyUtils</b>		CC0-1.0; Apache-2.0
<b>Log4j Implemented Over SLF4J</b>		Apache-2.0; MIT
<b>MapStruct Core</b>	Gunnar Morling, Filip Hrisafov	Apache-2.0
<b>mchange-commons-java</b>		LGPL-2.1-only; Multi-license: LGPL-2.1-only OR MIT
<b>metadata-extractor</b>		Apache-2.0
<b>Metrics Core</b>		CC0-1.0; Apache-2.0
<b>Netty/Buffer</b>		Apache-2.0
<b>Netty/Buffer</b>		Apache-2.0
<b>Netty/Codec</b>		Apache-2.0; Multi-license: Apache-2.0 OR cc-pd
<b>Netty/Codec</b>		Apache-2.0; Multi-license: Apache-2.0 OR cc-pd
<b>Netty/Codec/DNS</b>		Apache-2.0
<b>Netty/Codec/HAProxy</b>		Apache-2.0
<b>Netty/Codec/HTTP</b>		Apache-2.0; Multi-license: Apache-2.0 OR MIT; Multi-license: Apache-2.0 OR BSD-3-Clause
<b>Netty/Codec/HTTP2</b>		Apache-2.0
<b>Netty/Codec/Memcache</b>		Apache-2.0
<b>Netty/Codec/MQTT</b>		Apache-2.0
<b>Netty/Codec/Redis</b>		Apache-2.0
<b>Netty/Codec/SMTP</b>		Apache-2.0
<b>Netty/Codec/Socks</b>		Apache-2.0
<b>Netty/Codec/Stomp</b>		Apache-2.0
<b>Netty/Codec/XML</b>		Apache-2.0
<b>Netty/Common</b>		Apache-2.0; Multi-license: Apache-2.0 OR CC0-1.0; Multi-license: Apache-2.0 OR MIT
<b>Netty/Common</b>		Apache-2.0; Multi-license: Apache-2.0 OR CC0-1.0; Multi-license: Apache-2.0 OR MIT
<b>Netty/Handler</b>		Apache-2.0
<b>Netty/Handler</b>		Apache-2.0
<b>Netty/Handler/Proxy</b>		Apache-2.0

Netty/Resolver		Apache-2.0
Netty/Resolver		Apache-2.0
Netty/Resolver/DNS		Apache-2.0
Netty/Resolver/DNS/Classes/MacOS		Apache-2.0
Netty/Resolver/DNS/Native/MacOS		Apache-2.0
Netty/Transport		Apache-2.0
Netty/Transport		Apache-2.0
Netty/Transport/Classes/Epoll		Apache-2.0
Netty/Transport/Classes/KQueue		Apache-2.0
Netty/Transport/Native/Epoll		Apache-2.0
Netty/Transport/Native/Epoll		Apache-2.0
Netty/Transport/Native/KQueue		Apache-2.0
Netty/Transport/Native/Unix/Common		Apache-2.0
Netty/Transport/Native/Unix/Common		Apache-2.0
Netty/Transport/RXTX		Apache-2.0
Netty/Transport/SCTP		Apache-2.0
Netty/Transport/UDT		Apache-2.0
Nimbus Content Type		Apache-2.0
Nimbus JOSE+JWT		Apache-2.0
Nimbus LangTag		Apache-2.0
OAuth2.0 SDK with OpenID Connect extensions		Apache-2.0
Objenesis		Apache-2.0
OGNL - Object Graph Navigation Library		Apache-2.0; BSD-3-Clause
OGNL - Object Graph Navigation Library		Apache-2.0; BSD-3-Clause
Old JAXB Core		SAX-PD
Old JAXB Runtime		BSD-3-Clause; Multi-license: Apache-2.0 OR BSD-3-Clause OR MIT OR SAX-PD; Multi-license: BSD-3-Clause OR SAX-PD; SAX-PD
OXML schemas		Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR EPL-1.0 OR MIT; Apache-1.1
OpenSAML:: Core		Apache-2.0
OpenSAML:: Messaging API		Apache-2.0
OpenSAML:: Profile API		Apache-2.0
OpenSAML:: SAML Provider API		Apache-2.0
OpenSAML:: SAML Provider Implementations		Apache-2.0
OpenSAML:: Security API		Apache-2.0
OpenSAML:: Security Implementation		Apache-2.0
OpenSAML:: SOAP Provider API		Apache-2.0; W3C; W3C-19980720

OpenSAML::SOAP Provider Implementations		Apache-2.0
OpenSAML::Storage API		Apache-2.0
OpenSAML::XML Security API		Apache-2.0; W3C; W3C-19980720
OpenSAML::XML Security Implementation		Apache-2.0
org.apache.poi.xwpf.converter.core		MIT
org.opentest4j:opentest4j	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen	Apache-2.0
parso	Petr Tsurinov, Igor Printsev	Apache-2.0
PDFBox JBIG2 ImageIO plugin	Andreas Lehmkuhler, Adam Nichols, Ben Litchfield, Brian Carrier, Daniel Wilson, Guillaume Bailleul, Jörg Henne, Sebastian Holder, John Hewson, Jeremias Maerki, Kevin Jackson, Johannes Koch, Carolin Köhler, Matthäus Mayer, Maruan Sayhoun, Philipp Koch, Tim Allison, Thomas Chojecki, Timo Boehme, Tilman Hausherr	Apache-2.0
picocli		Apache-2.0
Prometheus Java Simpleclient		Apache-2.0; CC0-1.0
Prometheus Java Simpleclient Common		Apache-2.0
QDOX	Mike Royle, Peter Donald, James Strachan, Nick Pomfret, Chris Stevenson, Ben Hogan, Laurent Etienneble, Shawn Chain, Brian Slesinsky, James Lee, Eric Redmond, Robert Scholte, Joe Walnes, Aslak Hellesoy, Paul Hammant, Mike Williams, Mauro Talevi	Apache-2.0
SentimentAnalysisParser		Apache-2.0
Simple API for CSS		W3C-19980720
SLF4J API Module		MIT
SLF4J API Module		MIT
SLF4J API Module		MIT
SnakeYAML	Andrey Somov, Alexander Maslov, Jordan Angold	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR EPL-1.0 OR GPL-2.0-or-later OR LGPL-2.1-or-later
SnakeYAML	Andrey Somov, Alexander Maslov, Jordan Angold	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR EPL-1.0 OR GPL-2.0-or-later OR LGPL-2.1-or-later
SparseBitSet		Apache-2.0
Spring AOP		Apache-2.0
Spring AOP		Apache-2.0
Spring Beans		Apache-2.0
spring-boot		Apache-2.0
spring-boot-autoconfigure		Apache-2.0
spring-boot-starter		
spring-boot-starter-logging		
spring-boot-starter-tomcat		
Spring Commons Logging Bridge		Apache-2.0
Spring Context		Apache-2.0
Spring Core		Apache-2.0; BSD-3-Clause

Spring Expression Language (SpEL)		Apache-2.0
Spring Expression Language (SpEL)		Apache-2.0
springfox-schema	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-spi	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-spring-web	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-swagger-common	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
Spring JDBC		Apache-2.0
Spring JMS		Apache-2.0
spring-ldap-core	Eric Dalquist, Marius Scurtescu, Tim Terry, Keith Barlow, Paul Harvey, Marvin S. Addison, Rob Winch, Mattias Hellborg Arthursson, Ulrik Sandberg	Apache-2.0
Spring Plugin - Core		Apache-2.0
Spring Plugin - Metadata Extension		Apache-2.0
spring-security-crypto	Rob Winch, Joe Grandja	Apache-2.0; MIT
spring-security-oauth2-core	Rob Winch, Joe Grandja	Apache-2.0
Spring Transaction		Apache-2.0
Spring Web		Apache-2.0
Spring Web		Apache-2.0; public-domain
Spring Web MVC		Apache-2.0
Spring Web MVC		Apache-2.0
Stax2 API		Apache-2.0; ImageMagick; BSD-2-Clause
StAX API	Aleksander Slominski, Chris Fry	Apache-2.0
telemetry		Apache-2.0
telemetry-core		Apache-2.0
testng	Cedric Beust, Julien Herr, Krishnan Mahadevan	Apache-2.0
thymeleaf	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0; MIT
thymeleaf-layout-dialect	Justin Munn, Oliver Niekrenz	Apache-2.0
thymeleaf-spring4	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0; MIT
tomcat-annotations-api		Apache-2.0
tomcat-embed-core		Apache-2.0; Multi-license: CDDL-1.1 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; CDDL-1.1
tomcat-embed-el		Apache-2.0
tomcat-embed-websocket		Apache-2.0
TXW2 Runtime		Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception OR SAX-PD; Multi-license: CDDL-1.0 OR CDDL-1.1 OR GPL-2.0-with-classpath-exception; Apache-2.0
unbescape		Apache-2.0
unbescape		Apache-2.0
Vaadin Sass Compiler		Apache-2.0; Multi-license: Apache-2.0 OR W3C-19980720
vaadin-shared	Artur Signell, Marko Gronroos	Apache-2.0

<b>VT Crypt Library</b>	Daniel Fisher, Marvin Addison	Apache-2.0
<b>VT Dictionary Libraries</b>	Daniel Fisher, Marvin Addison	Apache-2.0
<b>Woodstox</b>		Apache-2.0; ImageMagick
<b>wstx-asl</b>		Apache-2.0; ImageMagick
<b>Xalan Java Serializer</b>		Apache-2.0; Multi-license: Apache-2.0 OR TCL; W3C-19980720
<b>Xerces2j</b>		
<b>XHTMLRenderer core-renderer</b>		LGPL-2.1-only; LGPL-2.1-or-later; GPL-2.0-or-later
<b>xml-apis</b>		Multi-license: W3C OR W3C-19980720; Apache-2.0; Multi-license: Apache-2.0 OR SAX-PD; SAX-PD; W3C-19980720; Multi-license: Apache-1.1 OR Apache-2.0; W3C
<b>XmlBeans</b>	Cezar Andrei, David Bau, Patrick Calahan, Yana Kadiyska, Ken Kress, Kevin Krouse, Laurence Moroney, Radu Preotiuc, Cliff Schmidt, Dutta Satadip, Eric Vasilik, David Waite, Scott Ziegler	Apache-2.0; Apache-1.1; Multi-license: Apache-2.0 OR IPA
<b>XML Commons External Components XML APIs Extensions</b>		Apache-2.0
<b>XMPCore Shaded</b>		Apache-2.0

[www.agilent.com/chem/agilentlims-licenses](http://www.agilent.com/chem/agilentlims-licenses)

This information is subject to change without notice.

5994-4865EN

© Agilent Technologies, Inc. 2022

Published in Europe, April 14, 2022

