

Agilent SLIMS 6.9

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 6.9.

Title	Authors	Declared License
Apache Ant Core		Apache-2.0; Multi-license: Apache-2.0 OR SAX-PD OR W3C; Info-ZIP
Apache Commons BeanUtils	Robert Burrell Donkin, Dion Gillard, Craig McClanahan, Geir Magnusson Jr., Scott Sanders, James Strachan, Rodney Waldhoff, Martin van den Bemt, Yoav Shapira, Niall Pemberton, Simon Kitching, James Carman, Benedikt Ritter, Tim O'Brien, David Eric Pugh, Rodney Waldhoff, Morgan James Delagrange, John E. Conlon, Stephen Colebourne, Gary Gregory	Apache-2.0
Apache Commons Codec	Henri Yandell, Tim O'Brien, Scott Sanders, Rodney Waldhoff, Daniel Rall, Jon S. Stevens, Gary Gregory, David Graham, Julius Davies, Thomas Neidhart, Rob Tompkins, Christopher O'Brien, Martin Redington, Jeffery Dever, Steve Zimmermann, Benjamin Walstrum, Oleg Kalnichevski, Dave Dribin, Alex Karasulu, Matthew Inger	Apache-2.0
Apache Commons Collections	Matt Benson, James Carman, Stephen Colebourne, Robert Burrell Donkin, Morgan Delagrange, Gary Gregory, Matthew Hawthorne, Dipanjan Laha, Geir Magnusson, Luc Maisonobe, Craig McClanahan, Thomas Neidhart, Adrian Nistor, Arun M. Thomas, Rodney Waldhoff, Henri Yandell, Rob Tompkins, Rafael U. C. Afonso, Max Rydahl Andersen, Avalon	Apache-2.0
Apache Commons Compress	Torsten Curdt, Stefan Bodewig, Sebastian Bazley, Christian Grobmeier, Julius Davies, Damjan Jovanovic, Emmanuel Bourg, Gary Gregory, Rob Tompkins, Peter Alfred Lee, Wolfgang Glas, Christian Kohlschütter, Bear Giles, Michael Kuss, Lasse Collin, John Kodis, BELUGA BEHR, Simon Spero, Michael Husegger, Arturo Bernal	Apache-2.0; Multi-license: Apache-2.0 OR Info-ZIP; bzip2-1.0.6
Apache Commons CSV	Henri Yandell, Martin van den Bemt, Yonik Seeley, Emmanuel Bourg, Gary Gregory, Benedikt Ritter, Rob Tompkins	Apache-2.0
Apache Commons DBCP	Todd Carmichael, Wayne Woodfield, Dain Sundstrom, Philippe Mouawad, Glenn L. Nielsen, James House, James Ring, Peter Wicks, Morgan Delagrange, 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
Apache Commons FileUpload	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, Dion Gillard	Apache-2.0
Apache Commons IO	Scott Sanders, Dion Gillard, Nicola Ken Barozzi, Henri Yandell, Stephen Colebourne, Jeremias Maerki, Matthew Hawthorne, Martin Cooper, Rob Oxspring, Jochen Wiedmann, Niall Pemberton, Jukka Zitting, Gary Gregory, Kristian Rosenvold, Rahul Akolkar, Jason Anderson, Nathan Beyer, Emmanuel Bourg, Chris Eldredge, Magnus Grimsell	Apache-2.0

Apache Commons IO	Scott Sanders, dlon Gillard, Nicola Ken Barozzi, Henri Yandell, Stephen Colebourne, Jeremias Maerki, Matthew Hawthorne, Martin Cooper, Rob Oxspring, Jochen Wiedmann, Niall Pemberton, Jukka Zitting, Gary Gregory, Kristian Rosenvold, Rahul Akolkar, Jason Anderson, Nathan Beyer, Emmanuel Bourg, Chris Eldredge, Magnus Grimsell	Apache-2.0
Apache 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, Morgan Delagrange, Ringo De Smet	Apache-2.0
Apache Commons Logging	Juozas Baliuka, Morgan Delagrange, Peter Donald, Robert Burrell Donkin, Simon Kitching, Dennis Lundberg, Costin Manolache, Craig McClanahan, Thomas Neidhart, Scott Sanders, Richard Sitze, Brian Stansberry, Rodney Waldhoff, Matthew P. Del Buono, Vince Eagen, Peter Lawrey, Berin Loritsch, Philippe Mouawad, Neeme Praks	Apache-2.0
Apache Commons Validator	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, Makoto Uchino	Apache-2.0
Apache FontBox		Apache-2.0; Multi-license: Apache-2.0 OR OFL-1.1; BSD-3-Clause
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrindle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0; Multi-license: Apache-2.0 OR public-domain; BSD-3-Clause; MIT
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrindle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache HttpClient		Apache-2.0
Apache HttpCore		Apache-2.0
Apache Log4j		Apache-2.0; Multi-license: Apache-2.0 OR TCL
Apache Log4j API		Apache-2.0; Multi-license: Apache-2.0 OR MIT
Apache Log4j Core		Apache-2.0; Multi-license: Apache-2.0 OR CDDL-1.0 OR MIT; Apache-1.1
Apache Log4j SLF4J Binding		Apache-2.0; Multi-license: Apache-2.0 OR MIT; Apache-1.1
Apache Log4j Spring Boot Support		Apache-2.0
Apache Log4j Web		Apache-2.0; Apache-1.1
Apache Lucene (module: core)		Apache-2.0; Multi-license: Apache-2.0 OR BSD-2-Clause OR BSD-3-Clause OR ICU OR MIT; BSD-3-Clause; Multi-license: MIT OR naist-2003 OR Unicode-DFS-2020; MIT; ICU
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: queryparser)		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
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		Unicode; Zlib; BSD-3-Clause; APAFML; OFL-1.1; Apache-2.0
Apache PDFBox		Unicode; Zlib; BSD-3-Clause; APAFML; OFL-1.1; Apache-2.0
Apache PDFBox Debugger		Apache-2.0; Apache-1.1
Apache PDFBox Debugger		Apache-2.0; Apache-1.1
Apache PDFBox tools		Apache-2.0; Apache-1.1
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 core		Apache-2.0; Apache-1.1
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 XML Security for Java	Sean Mullan, Berin Lautenbach, Axl Mattheus, Erwin van der Koog, Raul Benito Garcia, Vishal Mahajan, Milan Tomic, Colm O hEigearthaigh	Apache-2.0; Multi-license: Apache-2.0 OR MIT; W3C; W3C-19980720
Apache XML Security for Java	Sean Mullan, Berin Lautenbach, Axl Mattheus, Erwin van der Koog, Raul Benito Garcia, Vishal Mahajan, Milan Tomic, Colm O hEigearthaigh	Apache-2.0; Multi-license: Apache-2.0 OR MIT; W3C-19980720; Apache-1.1; W3C
ASM based accessors helper used by json-smart	Uriel Chemouni, ZhangJian He	Apache-2.0
AutoValue Annotations		Apache-2.0
barbecue	Sean C. Sullivan, Werner Keil	BSD-3-Clause
Barcode4J	Jeremias MÅrki, Web Maestro Clay Leeds, Nicola Ken Barozzi, Dietmar BÅrkle	Apache-2.0
batik-awt-util		Apache-2.0
batik-bridge		Apache-2.0
batik-dom		Apache-2.0
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 PKIX, CMS, EAC, TSP, PKCS, OCSP, CMP, and CRMF APIs	The Legion of the Bouncy Castle Inc.	BouncyCastle
Bouncy Castle Provider	The Legion of the Bouncy Castle Inc.	BouncyCastle; MIT
Bouncy Castle Provider	The Legion of the Bouncy Castle Inc.	BouncyCastle; MIT
Bouncy Castle S/MIME API	The Legion of the Bouncy Castle Inc.	BouncyCastle
Bouncy Castle S/MIME API	The Legion of the Bouncy Castle Inc.	BouncyCastle

c3p0		LGPL-2.1-only; Multi-license: LGPL-2.1-only OR EPL-1.0
Camel :: API		Apache-2.0
Camel :: Attachments		Apache-2.0
Camel :: Base Engine		Apache-2.0
Camel :: Core		Apache-2.0
Camel :: Core Model		Apache-2.0
Camel :: Core XML		Apache-2.0
Camel :: File		Apache-2.0
Camel :: HazelCast		Apache-2.0
Camel :: JMS		Apache-2.0
Camel :: Kafka		Apache-2.0
Camel :: Mail		Apache-2.0
Camel :: Main		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 :: Spring XML		Apache-2.0
Camel :: Support		Apache-2.0
ClassMate	Tatu Saloranta, Brian Langel	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	Daniel Rall, Stephen Colebourne, Henri Yandell, Steven Caswell, Robert Burrell Donkin, Gary D. Gregory, Phil Steitz, Fredrik Westermarck, James Carman, Niall Pemberton, Matt Benson, Joerg Schaible, Oliver Heger, Paul Benedict, C. Scott Ananian, Chris Audley, Stephane Bailliez, Michael Becke, Benjamin Bentmann, Ola Berg	Apache-2.0
Date Calculator Common		Apache-2.0
Date Calculator for Joda		Apache-2.0
dom4j		dom4j
domino-jackson		Apache-2.0
EventBinder		Apache-2.0
FindBugs-jsr305		Apache-2.0; CC-BY-2.5
fr.opensagres.poi.xwpf.converter.core		MIT
fr.opensagres.poi.xwpf.converter.pdf		MIT
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		Apache-2.0; Multi-license: Apache-2.0 OR CC0-1.0
Google Guice - Extensions - AssistedInject		Apache-1.1; Apache-2.0

Guava: Google Core Libraries for Java		Apache-2.0; CC0-1.0; Multi-license: CC0-1.0 OR cc-pd
Guava GWT compatible libs		Apache-2.0; Multi-license: Apache-2.0 OR CC0-1.0; CC0-1.0
gwt-servlet		Apache-2.0; CC0-1.0
gwt-user		Apache-2.0; public-domain
hazelcast		Multi-license: Apache-2.0 OR MIT; Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR CC0-1.0 OR CDDL-1.0 OR EPL-2.0 OR GPL-2.0-only OR JSON OR LGPL-2.0-only; JSON; MIT; Multi-license: BSD-3-Clause OR GPL-2.0-only OR MIT; GPL-3.0-only
hazelcast		Apache-2.0; MIT; AML; CC0-1.0
hazelcast-spring		Apache-2.0; BSD-3-Clause
hazelcast-spring		Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR GPL-2.0-only OR MIT; Multi-license: Apache-2.0 OR BSD-3-Clause OR CC0-1.0 OR CDDL-1.0 OR EPL-2.0 OR GPL-2.0-only OR JSON OR LGPL-2.0-only
hibernate-core - relocation	The Hibernate Development Team	LGPL-2.1-or-later; LGPL-3.0-only; EPL-2.0
hibernate-core - relocation	The Hibernate Development Team	EPL-2.0; LGPL-3.0-only; LGPL-2.1-or-later
Hibernate Validator Engine - Relocation Artifact		Apache-2.0
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 GPL-2.0-with-autoconf-exception OR ICU OR public-domain OR Unicode-DFS-2020; ICU; MIT; IPA
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 datatype: Guava	Steven Schlansker	Apache-2.0
Jackson datatype: jdk8		Apache-2.0
Jackson datatype: JSR310	Nick Williams	Apache-2.0
jakarta-regexp		Apache-2.0
JasperReports Font Extension	Teodor Danciu	LGPL-3.0-only
JasperReports Functions	Teodor Danciu	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
JavaBeans Activation Framework API jar		Multi-license: CDDL-1.0 OR GPL-2.0-only 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 GPL-2.0-with-classpath-exception; Apache-2.0
JavaBeans(TM) Activation Framework		Multi-license: CDDL-1.0 OR CDDL-1.1; Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; CDDL-1.0
java jwt	Auth0, Hernan Zalazar, Luciano Balmaceda	MIT
JavaMail API		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 cockroach 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
Java Servlet API	Ed Burns, Shing Wai Chan	Multi-license: CDDL-1.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Apache-2.0; Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: CDDL-1.1 OR GPL-2.0-with-classpath-exception
Javassist	Shigeru Chiba, Andrew Dinn, Kabir Khan, Scott Marlow	Multi-license: Apache-2.0 OR LGPL-2.1-only OR LGPL-2.1-or-later OR MPL-1.1 OR MPL-2.0; Multi-license: Apache-2.0 OR LGPL-2.1-or-later OR MPL-1.1
javax.annotation-api	Linda De Michiel	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
javax.inject		Apache-2.0
javax.persistence-api		Multi-license: BSD-3-Clause OR EPL-1.0; Apache-2.0
jaxb-api	Martin Grebac, Iaroslav Savitskiy	Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception OR GPL-3.0-only; Multi-license: 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; Apache-2.0
jaxen	Bob McWhirter, James Strachan, David Megginson, Erwin Bolwitt, Mark A. Belonga, Christian Nentwich, Alexander Day Chaffee, Jan Dvorak, Attila Szegedi, Peter Royal, Scott Sanders, Brian Ewins, Elliotte Rusty Harold, Ryan Gustafson, David Peterson, Mark Wilson, Jacob Krome, Michael Brennan, Jason Hunter, Brett McLaughlin	BSD-3-Clause; Apache-2.0
jaxen	Bob McWhirter, James Strachan, David Megginson, Erwin Bolwitt, Mark A. Belonga, Christian Nentwich, Alexander Day Chaffee, Jan Dvorak, Attila Szegedi, Peter Royal, Scott Sanders, Brian Ewins, Elliotte Rusty Harold, Ryan Gustafson, David Peterson, Mark Wilson, Jacob Krome, Michael Brennan, Jason Hunter, Brett McLaughlin	BSD-2-Clause; BSD-3-Clause
jboss-transaction-api_1.2_spec	Stephen Felts	Multi-license: GPL-2.0-with-classpath-exception OR public-domain
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
jCSV		Apache-2.0
JDOM	Jason Hunter, Rolf Lear	Multi-license: Apache-2.0 OR Saxpath; Saxpath
JGraphT - Core		Multi-license: EPL-1.0 OR EPL-2.0; Multi-license: EPL-1.0 OR EPL-2.0 OR MIT
Joda-Time	Stephen Colebourne, Brian S O'Neill, Sergi Almacellas Abellana, Adrian Acala, Afif Ahmed, AlexMe951, Guy Allard, Brad Arndt, Mikey Austin, Dominik Balazs, Oren Benjamin, Besnik Bleta, Fredrik Borgh, Dave Brosius, Dan Cavallaro, Luc Claes, Emiliano Claria, Dan Cojocar, Evgeniy Devyatikh, dspitfire	Apache-2.0; Multi-license: Apache-2.0 OR MIT
JSON.simple	Yidong Fang	Apache-2.0
JSON Small and Fast Parser	Uriel Chemouni, Eitan Raviv, ZhangJian He	Apache-2.0
jsoup Java HTML Parser	Jonathan Hedley	MIT

Kotlin Stdlib Jdk8	Kotlin Team	Apache-2.0
Kotlin Stdlib Jdk8	Kotlin Team	Apache-2.0
Kubernetes Discovery Plugin for Hazelcast	Christoph Engelbert (@noctarius2k)	Apache-2.0
Liquibase	Nathan Voxland	Apache-2.0
liquibase-slf4j	Matt Bertolini	MIT
Lucene Common Analyzers		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 IPA; Multi-license: Apache-2.0 OR BSD-3-Clause; BSD-3-Clause; Multi-license: MIT OR naist-2003 OR Unicode-DFS-2020
mariadb-java-client	mariadb jdbc developers	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
mariadb-java-client	mariadb jdbc developers	LGPL-2.1-only
mchange-commons-java		Multi-license: EPL-1.0 OR LGPL-2.1-only; LGPL-2.1-only; Multi-license: LGPL-2.1-only OR MIT
micrometer-core	Tommy Ludwig	Apache-2.0; Multi-license: Apache-2.0 OR MIT
micrometer-core	Tommy Ludwig	Apache-2.0; Multi-license: Apache-2.0 OR MIT
micrometer-registry-new-relic		Apache-2.0
micrometer-registry-prometheus	Tommy Ludwig	Apache-2.0
micrometer-registry-prometheus	Tommy Ludwig	Apache-2.0
Microsoft JDBC Driver for SQL Server		MIT; Apache-2.0
Microsoft JDBC Driver for SQL Server		MIT; Apache-2.0
Nimbus JOSE+JWT	Vladimir Dzhuvinov	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		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
OpenSAML :: Core		Apache-2.0
OpenSAML :: SAML Provider API		Apache-2.0
OpenSAML :: SAML Provider Implementations		Apache-2.0
OWASP Java HTML Sanitizer		BSD-3-Clause
quartz		Apache-2.0
RCaller	Mehmet Hakan Satman, M. Batchkarov, Paul Curcean, Kopilov Aleksandr	LGPL-3.0-only
SLF4J API Module		MIT
SnakeYAML	Andrey Somov, Alexander Maslov	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	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
Spring AOP	Juergen Hoeller	Apache-2.0
Spring Beans	Juergen Hoeller	Apache-2.0

spring-boot	Pivotal	Apache-2.0
spring-boot-devtools	Pivotal	Apache-2.0
spring-boot-starter	Pivotal	Apache-2.0
spring-boot-starter-log4j2	Pivotal	Apache-2.0
spring-boot-starter-thymeleaf	Pivotal	Apache-2.0
spring-boot-starter-tomcat	Pivotal	Apache-2.0
spring-boot-starter-web	Pivotal	Apache-2.0
Spring Commons Logging Bridge	Juergen Hoeller	Apache-2.0
Spring Context	Juergen Hoeller	Apache-2.0
Spring Context Support	Juergen Hoeller	Apache-2.0
Spring Core	Juergen Hoeller	Apache-2.0
springfox-core	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-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 JDBC	Juergen Hoeller	Apache-2.0
Spring JMS	Juergen Hoeller	Apache-2.0
spring-ldap-core	Pivotal	Apache-2.0
Spring Messaging	Juergen Hoeller	Apache-2.0
Spring Object/Relational Mapping	Juergen Hoeller	Apache-2.0
spring-security-config	Pivotal	Apache-2.0
spring-security-core	Pivotal	Apache-2.0
spring-security-crypto	Pivotal	Apache-2.0; MIT
spring-security-ldap	Pivotal	Apache-2.0
spring-security-oauth2-authorization-server	Joe Grandja, Steve Riesenberg	Apache-2.0
spring-security-oauth2-client	Pivotal	Apache-2.0
spring-security-oauth2-core	Pivotal	Apache-2.0
spring-security-oauth2-jose	Pivotal	Apache-2.0
spring-security-oauth2-resource-server	Pivotal	Apache-2.0
spring-security-saml2-service-provider	Pivotal	Apache-2.0
spring-security-web	Pivotal	Apache-2.0
Spring Transaction	Juergen Hoeller	Apache-2.0
Spring Web	Juergen Hoeller	Apache-2.0
Spring Web MVC	Juergen Hoeller	Apache-2.0
Stax2 API		Apache-2.0; ImageMagick
StAX API	Aleksander Slominski, Chris Fry	Apache-2.0
swagger-annotations		Apache-2.0
swagger-models		Apache-2.0
thymeleaf	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0
thymeleaf-spring5	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0
tomcat-annotations-api		Apache-2.0
tomcat-embed-core		Apache-2.0; Multi-license: Apache-2.0 OR CDDL-1.0; Multi-license: CDDL-1.1

		OR GPL-2.0-with-classpath-exception; CC0-1.0; Multi-license: CDDL-1.0 OR GPL-2.0-with-classpath-exception
tomcat-embed-jasper		Apache-2.0
vaadin-server	Artur Signell, Marko Gronroos	Apache-2.0; atmosphere-0.4; Multi-license: Apache-2.0 OR atmosphere-0.4
vaadin-shared	Artur Signell, Marko Gronroos	Apache-2.0; Multi-license: Apache-2.0 OR atmosphere-0.4
vaadin-spring		Apache-2.0
vaadin-themes	Artur Signell, Marko Gronroos	Apache-2.0
VT Password Libraries	Daniel Fisher, Marvin Addison	Apache-2.0
Woodstox	Tatu Saloranta	Apache-2.0; ImageMagick
Xalan Java		Apache-2.0; W3C-19980720; Multi-license: Apache-2.0 OR TCL
Xerces2-j	Apache Software Foundation	Apache-2.0; Multi-license: Apache-2.0 OR IPA; W3C-19980720
Xerces2-j	Apache Software Foundation	W3C-19980720; Apache-2.0; Multi-license: Apache-2.0 OR IPA
xml-apis		Multi-license: Apache-2.0 OR SAX-PD OR unknown; Apache-2.0; Multi-license: W3C OR W3C-19980720; SAX-PD; Multi-license: Apache-2.0 OR SAX-PD; W3C-19980720; W3C
ZXing Core		Apache-2.0
Deep Dependencies		
Title	Authors	Declared License
Adobe XMPCore		BSD-2-Clause
AntLR		BSD-2-Clause
AOP alliance		public-domain
Apache Ant + ANTLR		Apache-2.0; Multi-license: Apache-2.0 OR SAX-PD OR W3C; Apache-1.1
Apache Ant + JUnit		Apache-2.0; Multi-license: Apache-2.0 OR SAX-PD OR W3C; Apache-1.1
Apache Ant Launcher		Apache-2.0; Multi-license: Apache-2.0 OR SAX-PD OR W3C; Apache-1.1
Apache Commons CLI	James Strachan, Bob McWhirter, John Keyes, Rob Oxspring, Emmanuel Bourg, Thomas Neidhart, Rob Tompkins, Gary Gregory, Beluga Behr, Peter Donald, Brian Egge, Duncan Jones, Berin Loritsch, Peter Maddocks, Alexandru Mocanu, Andrew Shirley, Greg Thomas, Slawek Zachcial, Rubin Simons	Apache-2.0
Apache Commons CLI	James Strachan, Bob McWhirter, John Keyes, Rob Oxspring, Emmanuel Bourg, Thomas Neidhart, Rob Tompkins, Gary Gregory, Beluga Behr, Peter Donald, Brian Egge, Duncan Jones, Berin Loritsch, Peter Maddocks, Alexandru Mocanu, Andrew Shirley, Greg Thomas, Slawek Zachcial, Rubin Simons	Apache-2.0
Apache Commons Collections	Matt Benson, James Carman, Stephen Colebourne, Robert Burrell Donkin, Morgan Delagrang, Gary Gregory, Matthew Hawthorne, Dipanjan Laha, Geir Magnusson, Luc Maisonobe, Craig McClanahan, Thomas Neidhart, Adrian Nistor, Arun M. Thomas, Rodney Waldhoff, Henri Yandell, Rob Tompkins, Rafael U. C. Afonso, Max Rydahl Andersen, Avalon	Apache-2.0
Apache Commons Compress	Torsten Curdt, Stefan Bodewig, Sebastian Bazley, Christian Grobmeier, Julius Davies, Damjan Jovanovic, Emmanuel Bourg, Gary Gregory, Rob Tompkins, Peter Alfred Lee, Wolfgang Glas, Christian Kohlschütter, Bear	Apache-2.0; Multi-license: Apache-2.0 OR Info-ZIP

	Giles, Michael Kuss, Lasse Collin, John Kodis, BELUGA BEHR, Simon Spero, Michael Hausegger, Arturo Bernal	
Apache Commons Exec	Niklas Gustavsson, Benjamin Bentmann, Marco Ferrante, Jerome Lacoste, Milos Kleint, Pablo Hoertner, Niall Pemberton, Brett Porter, Trygve Laugst�, Siegfried Goeschl, Sebastian Bazley, Gary Gregory	Apache-2.0
Apache Commons Logging	Juozas Baliuka, Morgan Delagrance, Peter Donald, Robert Burrell Donkin, Simon Kitching, Dennis Lundberg, Costin Manolache, Craig McClanahan, Thomas Neidhart, Scott Sanders, Richard Sitze, Brian Stansberry, Rodney Waldhoff, Matthew P. Del Buono, Vince Eagen, Peter Lawrey, Berin Loritsch, Philippe Mouawad, Neeme Praks	Apache-2.0
Apache Commons Math	Mikkel Meyer Andersen, Bill Barker, S�@bastien Brisard, Albert Davidson Chou, Mark Diggory, Robert Burrell Donkin, Otmar Ertl, Luc Maisonobe, Tim O'Brien, J. Pietschmann, Dimitri Pourbaix, Gilles Sadowski, Greg Sterijevski, Brent Worden, Thomas Neidhart, Evan Ward, Eldar Agalarov, Tim Allison, C. Scott Ananian, Mark Anderson	Apache-2.0; BSD-3-Clause; BSD-2-Clause; Multi-license: Apache-2.0 OR BSD-3-Clause
Apache Commons Pool	Morgan Delagrance, Geir Magnusson, Craig McClanahan, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Robert Burrell Donkin, Sandy McArthur, Simone Tripodi, Gary Gregory, Matt Sicker	Apache-2.0; Multi-license: Apache-2.0 OR cc-pd; Apache-1.1
Apache FontBox		Apache-2.0; BSD-3-Clause; Apache-1.1
Apache Geronimo JMS Spec 2.0		Apache-2.0; Apache-1.1
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrindle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrindle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrindle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrindle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrindle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrindle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrindle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0; Multi-license: Apache-2.0 OR public-domain

	McCrinkle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrinkle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrinkle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrinkle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrinkle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrinkle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause
Apache Groovy	Guillaume Laforge, bob mcwhirter, James Strachan, Joe Walnes, Chris Stevenson, Jamie McCrinkle, Matt Foemmel, Alex Tkachman, Roshan Dawrani, Sam Pullara, Kasper Nielsen, Travis Kay, Zohar Melamed, John Wilson, Chris Poirier, Christiaan ten Klooster, Steve Goetze, Bing Ran, Jeremy Rayner, John Stump	Apache-2.0
Apache HttpClient Mime		Apache-2.0
Apache HttpClientCore		Apache-2.0
Apache JempBox		Apache-2.0
Apache Kafka		Apache-2.0; Multi-license: BSD-3-Clause OR EPL-2.0; BSD-3-Clause
Apache Log4j JUL Adapter		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 Velocity - Engine		Multi-license: Apache-2.0 OR MIT; Apache-2.0
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
ASM Core	Eric Bruneton, Eugene Kuleshov	BSD-3-Clause
attoparser	Daniel Fernandez	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
batik-shared-resources		Apache-2.0
batik-svg-dom		Apache-2.0
Bean Validation API	Emmanuel Bernard, Hardy Ferentschik, Gunnar Morling, Guillaume Smet	Apache-2.0
Bouncy Castle ASN.1 Extension and Utility APIs	The Legion of the Bouncy Castle Inc.	BouncyCastle
Bouncy Castle PKIX, CMS, EAC, TSP, PKCS, OCSP, CMP, and CRMF APIs	The Legion of the Bouncy Castle Inc.	BouncyCastle
Byte Buddy (without dependencies)		Apache-2.0
c3p0		LGPL-2.1-only; Multi-license: LGPL-2.1-only OR EPL-1.0
Camel :: Base		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 Processor		Apache-2.0
Camel :: Core Reifier		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 :: Health		Apache-2.0
Camel :: Language		Apache-2.0
Camel :: Log		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 :: 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
CGLIB	Juozas Baliuka, Chris Nokleberg, Matas Ramoska, Sam Berlin, Wes Biggs	Apache-2.0; BSD-3-Clause
Checker Qual	Michael Ernst, Suzanne Millstein	MIT
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 Pool	Morgan Delagrang, Geir Magnusson, Craig McClanahan, Rodney Waldhoff, David Weinrich, Dirk Verbeeck, Robert Burrell Donkin, Sandy McArthur, Gary Gregory, Todd Carmichael	Apache-2.0
Cryptacular Library	Daniel Fisher, Marvin S. Addison	Multi-license: Apache-2.0 OR LGPL-3.0-only
curvesapi	stormdollar, Dustin Spicuzza	Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause; Apache-2.0
dec		MIT
domino-jackson-super		Apache-2.0
Eclipse Compiler for Java(TM)		Multi-license: EPL-1.0 OR EPL-2.0; EPL-2.0; EPL-1.0
Elemental2 Core	J2CL Team	Apache-2.0
Elemental2 Dom	J2CL Team	Apache-2.0
Elemental2 Promise		Apache-2.0
error-prone annotations		Apache-2.0
Flute		W3C; W3C-19980720
fr.opensagres.poi.xwpf.converter.xhtml		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 InternalFutureFailureAccess and InternalFutures		Apache-2.0
Guava ListenableFuture only		Apache-2.0
GWT CLDR	The GWT Project Authors	Apache-2.0
GWT DateTimeFormat		Apache-2.0

gwt-elemental		Apache-2.0
GWT i18n		Apache-2.0
Hamcrest	Joe Walnes, Nat Pryce, Steve Freeman	Multi-license: BSD-2-Clause OR BSD-3-Clause; BSD-3-Clause
Hamcrest Core		BSD-3-Clause-Clear
HdrHistogram	Gil Tene	Multi-license: BSD-2-Clause OR CC0-1.0 OR public-domain; CC0-1.0; Multi-license: BSD-2-Clause OR CC0-1.0
Hibernate Commons Annotations	The Hibernate Development Team	LGPL-2.1-or-later
ISO Parser		Apache-2.0; EPL-1.0
istack common utility code runtime		Multi-license: BSD-3-Clause OR EPL-1.0 OR GPL-2.0-with-classpath-exception; BSD-3-Clause; Multi-license: Apache-2.0 OR BSD-3-Clause OR MIT OR SAX-PD; Multi-license: EPL-2.0 OR GPL-2.0-with-classpath-exception
J2ObjC Annotations		Apache-2.0
Jackson BOM	Tatu Saloranta	Apache-2.0
jackson-databind		Apache-2.0
Jackson-module-parameter-names		Apache-2.0
Jakarta Activation		Multi-license: BSD-3-Clause OR EPL-1.0 OR GPL-2.0-with-classpath-exception OR unknown; BSD-3-Clause; Multi-license: EPL-2.0 OR GPL-2.0-with-classpath-exception
Jakarta Activation API		Multi-license: BSD-3-Clause OR EPL-1.0 OR GPL-2.0-with-classpath-exception OR unknown; BSD-3-Clause; Multi-license: EPL-2.0 OR GPL-2.0-with-classpath-exception
Jakarta Annotations API	Linda De Michiel, Dmitry Kornilov	Multi-license: EPL-1.0 OR EPL-2.0 OR GPL-2.0-with-classpath-exception OR GPL-3.0-only; Multi-license: EPL-2.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception; Multi-license: BSD-3-Clause OR EPL-2.0 OR GPL-2.0-with-classpath-exception; Multi-license: EPL-2.0 OR GPL-2.0-with-classpath-exception
Jakarta Mail API	Jakarta XML Binding API	BSD-3-Clause; Multi-license: EPL-2.0 OR GPL-2.0-only OR GPL-2.0-with-classpath-exception
		Multi-license: BSD-3-Clause OR EPL-1.0 OR GPL-2.0-with-classpath-exception; 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
jandex		Apache-2.0
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
JavaBeans Activation Framework		Multi-license: CDDL-1.0 OR GPL-2.0-only 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 GPL-2.0-with-classpath-exception; Apache-2.0
JavaBeans(TM) Activation Framework		CDDL-1.0
Java Native Access	Timothy Wall, Matthias Blasing	Multi-license: Apache-2.0 OR LGPL-2.1-or-later
javaparser-core		LGPL-3.0-only; Multi-license: Apache-2.0 OR LGPL-3.0-only; LGPL-3.0-or-later;

		Multi-license: Apache-2.0 OR LGPL-3.0-or-later
java-support		Apache-2.0
JAXB Runtime		SPL-1.0; GPL-2.0-with-classpath-exception
JBoss Logging 3		Multi-license: Apache-2.0 OR public-domain; Apache-2.0
JCIP Annotations under Apache License		Apache-2.0
JCL 1.2 implemented over SLF4J		Apache-2.0
jcommander	Cedric Beust	Apache-2.0
JetBrains Java Annotations	JetBrains Team	Apache-2.0
JFreeChart		LGPL-3.0-only; LGPL-2.1-or-later; BSD-3-Clause
JHeaps		Apache-2.0
JLine	Guillaume Nodet	BSD-3-Clause; Apache-2.0
JMatIO		BSD-2-Clause
Joda-Time	Stephen Colebourne, Brian S O'Neill, Sergi Almacellas Abellana, Adrian Acala, Afif Ahmed, AlexMe951, Guy Allard, Brad Arndt, Mikey Austin, Dominik Balazs, Oren Benjamin, Besnik Blea, Fredrik Borgh, Dave Brosius, Dan Cavallaro, Luc Claes, Emiliano Claria, Dan Cojocar, Evgeniy Devyatykh, dspitfire	Apache-2.0
Joda-Time	Stephen Colebourne, Brian S O'Neill, Sergi Almacellas Abellana, Adrian Acala, Afif Ahmed, AlexMe951, Guy Allard, Brad Arndt, Mikey Austin, Dominik Balazs, Oren Benjamin, Besnik Blea, Fredrik Borgh, Dave Brosius, Dan Cavallaro, Luc Claes, Emiliano Claria, Dan Cojocar, Evgeniy Devyatykh, dspitfire	Apache-2.0; X11
jquery	James Ward	MIT
JsInterop Annotations		Apache-2.0
JsInterop Base		Apache-2.0
JTidy		W3C-19980720; MIT
JUL to SLF4J bridge		MIT
JUnit	David Saff, Kevin Cooney, Stefan Birkner, Marc Philipp, JUnit contributors	EPL-1.0
JUnit 5 (Bill of Materials)	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0
JUnit Jupiter API	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0
JUnit Jupiter Engine	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0
JUnit Platform Commons	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0
JUnit Platform Engine API	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0
JUnit Platform Launcher	Stefan Bechtold, Johannes Link, Marc Philipp, Matthias Merdes, Sam Brannen, Christian Stein, Juliette de Rancourt	EPL-1.0; EPL-2.0
Kotlin Stdlib	Kotlin Team	Apache-2.0
Kotlin Stdlib Common	Kotlin Team	Apache-2.0
Kotlin Stdlib Jdk7	Kotlin Team	Apache-2.0
LatencyUtils		Multi-license: CC0-1.0 OR public-domain; CC0-1.0; Apache-2.0

LZ4 and xxHash	Adrien Grand, Rei Odaïra	Apache-2.0
MapStruct Core	Gunnar Morling, Filip Hrisafov	Apache-2.0
mchange-commons-java		LGPL-2.1-only; Multi-license: LGPL-2.1-only OR MIT
metadata-extractor	Drew Noakes	Apache-2.0
Metrics Core		Apache-2.0
Netty/Buffer		Apache-2.0
Netty/Codec		Apache-2.0
Netty/Common		Apache-2.0
Netty/Handler		Apache-2.0
Netty/Resolver		Apache-2.0
Netty/Transport		Apache-2.0
Netty/Transport/Classes/Epoll		Apache-2.0
Netty/Transport/Native/Epoll		Apache-2.0
Netty/Transport/Native/Unix/Common		Apache-2.0
Nimbus Content Type		Apache-2.0
Nimbus LangTag	Vladimir Dzhuvinov	Apache-2.0
OAuth 2.0 SDK with OpenID Connect extensions	Vladimir Dzhuvinov	Apache-2.0
OGNL - Object Graph Navigation Library		Apache-2.0; BSD-3-Clause
Old JAXB Core		SAX-PD
Old JAXB Runtime		Multi-license: EPL-1.0 OR GPL-2.0-with-classpath-exception; Multi-license: Apache-2.0 OR BSD-3-Clause OR MIT OR SAX-PD; BSD-3-Clause; Multi-license: BSD-3-Clause OR SAX-PD; SAX-PD; Multi-license: EPL-2.0 OR GPL-2.0-with-classpath-exception
OOXML schemas		Apache-2.0; Multi-license: Apache-2.0 OR BSD-3-Clause OR EPL-1.0 OR MIT; Apache-1.1
OpenSAML :: Messaging API		Apache-2.0
OpenSAML :: Profile API		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.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, Jürg Henne, Sebastian Holder, John Hewson, Jeremias Maerki, Kevin Jackson, Johannes Koch, Carolin Köhler, Matthias Mayer, Maruan Sayhoun, Phillipp Koch, Tim Allison, Thomas Chojecki, Timo Boehme, Tilman Hausherr	Apache-2.0
picocli	Remko Popma	Apache-2.0
Prometheus Java Simpleclient	Brian Brazil	Apache-2.0; CC0-1.0

Prometheus Java Simpleclient Common	Brian Brazil	Apache-2.0
Prometheus Java Span Context Supplier - Common		Apache-2.0
Prometheus Java Span Context Supplier - OpenTelemetry		Apache-2.0
Prometheus Java Span Context Supplier - OpenTelemetry Agent		Apache-2.0
QDox	Robert Scholte, Joe Walnes, Aslak Hellesoy, Paul Hammant, Mike Williams, Mauro Talevi, Mike Royle, Peter Donald, James Strachan, Nick Pomfret, Chris Stevenson, Ben Hogan, Laurent Etiemble, Shawn Chain, Brian Slesinsky, James Lee, Eric Redmond	Apache-2.0
SentimentAnalysisParser		Apache-2.0
Simple API for CSS		W3C-19980720
snappy-java	Taro L. Saito	Apache-2.0
SparseBitSet		Apache-2.0
spring-boot-autoconfigure	Pivotal	Apache-2.0
spring-boot-starter-json	Pivotal	Apache-2.0
Spring Expression Language (SpEL)	Juergen Hoeller	Apache-2.0
springfox-schema	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
springfox-swagger-common	Marty Pitt, Dilip Krishnan, Adrian Kelly	Apache-2.0
spring-ldap-core	Pivotal	Apache-2.0
Spring Plugin - Core	Oliver Drotbohm	Apache-2.0
Spring Plugin - Metadata Extension		Apache-2.0
Stax2 API		Apache-2.0; ImageMagick; BSD-2-Clause
telemetry		Apache-2.0
telemetry-core	New Relic	Apache-2.0
testng	Cedric Beust, Julien Herr, Krishnan Mahadevan	Apache-2.0
thymeleaf-extras-java8time	Daniel Fernandez, Jose Miguel Samper, Emanuel Rabina	Apache-2.0
tomcat-embed-core		Apache-2.0
tomcat-embed-el		Apache-2.0
tomcat-embed-websocket		Apache-2.0
TXW2 Runtime		GPL-2.0-with-classpath-exception; SPL-1.0
unbescape		Apache-2.0
Vaadin Sass Compiler		Apache-2.0; Multi-license: Apache-2.0 OR W3C-19980720
VT Crypt Library	Daniel Fisher, Marvin Addison	Apache-2.0
VT Dictionary Libraries	Daniel Fisher, Marvin Addison	Apache-2.0
Woodstox	Tatu Saloranta	Apache-2.0; ImageMagick
Xalan Java Serializer		Apache-2.0; Multi-license: Apache-2.0 OR TCL; W3C-19980720
XHTMLRenderer core-renderer		LGPL-2.1-only; LGPL-2.1-or-later; GPL-2.0-or-later
XmlBeans	POI Team	Apache-2.0; Multi-license: Apache-2.0 OR IPA; Apache-1.1
XML Commons External Components XML APIs Extensions		Apache-2.0
XMPCore Shaded		Apache-2.0
zstd-jni	Luben Karavelov	BSD-2-Clause

Learn more:

www.agilent.com/chem/agilentlims

U.S. and Canada

1-800-227-9770 agilent_inquiries@agilent.com

Europe

info_agilent@agilent.com

Asia Pacific

inquiry_lsca@agilent.com

This information is subject to change without notice.

DE14665644

© Agilent Technologies, Inc. 2022
Published in Europe, December 7, 2022
5994-5587EN

