

CONSISTENT QUALITY, MAXIMUM PRODUCTIVITY

Why gamble with your results? Agilent vials are the *only* vials that deliver time-saving, and cost-saving, advantages like these:

30+
Inspection points. So you get the tightest dimensional specifications, every time

10+
Instrument brands with which Agilent vials/caps are compatibility tested

127 (and growing)
Countries we deliver to across the world, from Albania to Vietnam

100s of millions
Agilent vials shipped worldwide every year

50%
Faster crimp speed: Our electric crimper lets you crimp your vial, not your style

33/51 Best in glass: All vials are made of type 33-51 coefficient of expansion for top performance

30%
Time savings using our full range of short thread screw top vials and caps

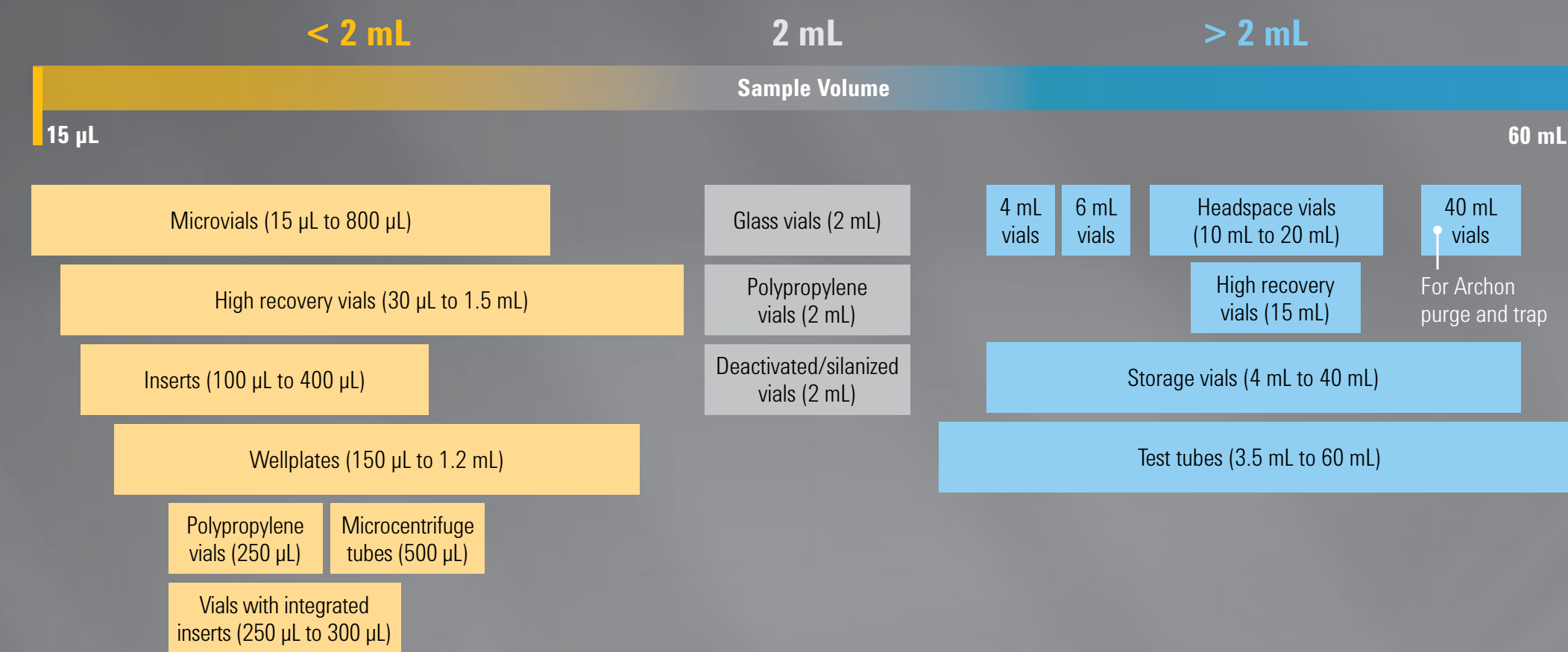
120 meters
Vials are small... but are manufactured in a facility as big as an aircraft hangar!

For an in-depth look at the Agilent Vials portfolio, including product brochure, crimping video, and white papers, visit www.agilent.com/chem/vialsresources

Information, descriptions, and specifications in this publication are subject to change without notice.
© Agilent Technologies, Inc. 2016
Printed in the U.S.A. June 16, 2016
5991-650EN

The industry's largest selection of sample containment products

The optimal sample size can be a function of many things, including analysis type, analytical platform, and sample availability. Agilent vials offer the same consistent performance across the entire size range, from 15 µL to 60 mL. What's more, they are manufactured to perform seamlessly with a variety of analytical instruments—regardless of make or model.



A grid of vial images with their dimensions and part numbers. The grid is organized into several sections:

- Inserts for Narrow Opening Vials (8 mm)**: 28 mm x 4.8 mm (150 µL, 5183-2088), 31 mm x 4.8 mm (150 µL, 5183-2089), 31 mm x 4.8 mm (200 µL, 5183-2090)
- Inserts for Wide Opening Vials (11 mm & 9 mm)**: 30 mm x 5.6 mm (250 µL, 5181-1270), 31 mm x 5.6 mm (250 µL, 5183-2085), 31 mm x 5.6 mm (400 µL, 5181-3377)
- Polypropylene Vials**: 32 mm x 12 mm (250 µL, 5188-2788), 32 mm x 12 mm (250 µL, 9301-0977 Glass Insert), 32 mm x 12 mm (700 µL, 5182-0567), 32 mm x 12 mm (250 µL, 5190-2242), 32 mm x 12 mm (250 µL, 5188-5390 Glass Insert)
- Narrow Opening Screw Top Vials (9 mm)**: 32 mm x 12 mm (250 µL, 5188-6591), 32 mm x 12 mm (1.2 mL, 5183-2030), 32 mm x 12 mm (1.3 mL, 5184-3550), 32 mm x 12 mm (1.5 mL, 5182-0714)
- Wide Opening Screw Top Vials (8 mm)**: 32 mm x 12 mm (1.5 mL, 5183-4428)
- Wide Opening Crimp Top Vials (11 mm)**: 32 mm x 12 mm (250 µL, 9301-1388), 32 mm x 12 mm (1.2 mL, 5182-3454), 32 mm x 12 mm (1.3 mL, 5184-3551), 32 mm x 12 mm (1.5 mL, 5181-3375)
- Large Volume Vials**: 45 mm x 15 mm (4 mL, 5183-4448), 37 mm x 22 mm (8 mL, 9301-1377), 37 mm x 22 mm (6 mL, 9301-1419)
- Headspace Vials**: 45 mm x 22 mm (10 mL, 5183-4475), 45 mm x 22 mm (20 mL, 5182-0838), 75 mm x 22 mm (20 mL, 5183-4474), 75 mm x 22 mm (20 mL, 5182-0837)

This represents a selection of vials. Please review the full portfolio on our website or see the new vials brochure, which contains the complete amber and clear selection.

Choose the right closure for your sample

Always make sure the septa you select are chemically compatible with your sample and solvent. Use this chart as a guide, but remember that chemical compatibility can vary based on solvent concentration, molecular weight, and temperature.

Septa Chemical Compatibility	PTFE	PTFE/Silicone	PTFE/Silicone/PTFE*	PTFE/Red Rubber	PTFE/Butyl
Acetonitrile	•	•	•	•	•
Hydrocarbons (hexane, heptane, methane)	•	•	•	•	•
Methanol	•	•	•	•	•
Benzene	•	•	•	•	•
THF	•	•	•	•	•
Toluene	•	•	•	•	•
DMF	•	•	•	•	•
DMSO	•	•	•	•	•
Ether	•	•	•	•	•
Chlorinated Solvents (methylene chloride)	•	•	•	•	•
Alcohols (ethanol)	•	•	•	•	•
Acetic Acid	•	•	•	•	•
Acetone	•	•	•	•	•
Phenol	•	•	•	•	•
Cyclohexane	•	•	•	•	•

*PTFE/silicone/PTFE has the same chemical compatibility of PTFE ONLY UNTIL PUNCTURED.

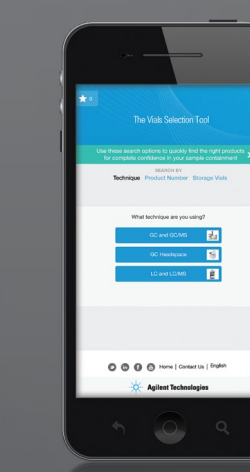
Use this chart to determine the right cap and septa combination, based on your application. Note: septa that are too thick can prevent the cap from fitting properly on the vial.

Cap and Septa Compatibility	High Performance Septa	Thin PTFE	PTFE/Silicone*	PTFE/Silicone/PTFE*	PTFE/Red Rubber	PTFE/Butyl
Part number ¹	5190-3986 (18 mm) 5190-3987 (20 mm)	5062-3582 (11 mm)	5190-7021 (9 mm)** 5190-7023 (9 mm pre-slit)**	5182-0723 (9 mm)	5181-1210 (11 mm)	5183-4479 (20 mm)
Temperature range	40 to 300 °C for up to 1 hour	Up to 260 °C	-40 °C to 200 °C	-40 °C to 200 °C	-40 °C to 90 °C	-50 °C to 150 °C
Use for multiple injections	No	No	Yes	Yes	No	No
Price	Most expensive	Very economical	Economical	Most expensive	Very economical	Economical
Resistance to coring	Excellent	None	Excellent	Excellent	None	None
Recommended for storage	No	No	Yes	Yes	No	No
Best for	High temperature headspace applications	Superior chemical inertness, short cycle times, and single injections	Most common HPLC and GC analyses, not as resistant to coring as P/S/P	Superior performance for ultra analysis, repeat injections, internal standards	Chlorosilanes more economical option for single injections	Organic solvents, acetic acids, impermeable to gasses

* Agilent silicone is peroxide cured, making it more inert and less likely to interact with samples.

** Now available in bonded.

¹ These are just a few options, many more are available.



The right vial is only a few clicks away

Use our online selection tool to quickly find the right products for complete confidence in your sample containment.

- Answer a few simple questions to identify your best options
- Search by technique, product number, vial type, or instrument manufacturer
- Make a perfect choice from more than 600 vials, caps, and septa

Go to www.agilent.com/chem/selectvials

Agilent CrossLab

From Insight to Outcome

Agilent CrossLab, the world leader in innovative laboratory services, software, and consumables, delivers vital, actionable insights to drive improved economic, operational, and scientific outcomes. Visit www.agilent.com/chem/crosslab