

Agilent QuickProbe GC/MS Supplies



Agilent supplies for Agilent instruments

Agilent is committed to optimizing your laboratory's productivity. We have produced this list of the most commonly ordered supplies and parts for the QuickProbe. Keep this list handy so you can quickly and easily find the supplies you need and minimize instrument downtime.

Insertion Probes	Part Number
QuickProbe probe holder For use with round tip and pocket tip probes	G3971-60200
Insertion probe round tip (100/pk) Recommended for liquids, dissolved samples, and tablets	5190-5118
Insertion probe pocket tip (100/pk) Recommended for powders	5190-5113
TSP probe holder For use with TSP solid probe vials	G3971-20251
TSP solid probe vials (100/pk) Recommended for solids	5190-3187
Additional Consumables	Part Number
External split vent cartridge kit Includes trap and 3 cartridges	RDT-1020
External split vent replacement cartridges (3/pk)	RDT-1023
Gas Clean carrier gas kit Includes one 1-position 1/8 in connecting unit (part number CP7988) and two carrier gas purifiers (part number CP17973)	CP17976
Gas Clean carrier gas purifier, 1/pk	CP17973

QuickProbe Inlet Consumables	Part Number
QuickProbe liner, UI, 4 mm id, frit	5190-5104
GC inlet seal, gold plated, with washer	5188-5367
QuickProbe Columns and Ferrules	Part Number
Agilent J&W DB-1ht QuickProbe GC column, 2 m, 0.25 mm, 0.10 μm, for use in QuickProbe Fast GC; high-temperature columns are highly recommended with the QuickProbe*	G3903-61006
Agilent J&W DB-1ms Ultra Inert QuickProbe GC column, 1 m, 0.18 mm, 0.18 μm , for use when connecting GC to MS	G3903-61007
Flexible Metal ferrule, UltiMetal Plus, 0.4 mm id, for 0.1 to 0.25 mm id fused silica tubing, 10/pk	G3188-27501
Ferrule, 0.4 mm id, 15% graphite/85% Vespel, 0.1 to 0.25 mm column, 10/pk	5181-3323
Ultimate Union kit	G3182-61580
Internal nut, CFT capillary fitting	G2855-20530
Self Tightening column nut, for Agilent mass spec interface transfer line	5190-5233

*Additional QuickProbe column phases can be ordered via the custom column process. www.agilent.com/chem/gccustomcolumns





Agilent supplies-to help you get the results you need

For greater confidence in your chromatography results, you can optimize your total system with quality Agilent columns and supplies. Our columns, supplies, kits, and accessories are designed, manufactured, and tested to rigorous Agilent specifications, under a quality system registered to ISO 9001. Why risk compromising your analytical results with anything less than genuine Agilent supplies?

When choosing your columns, Agilent J&W GC columns offer the lowest bleed levels, the best inertness for acids/bases/mixed functional compounds, and the tightest column-to-column reproducibility.

For a complete listing of Agilent columns and supplies, visit www.agilent.com/chem/supplies

www.agilent.com/chem/quickprobe

This information is subject to change without notice.

© Agilent Technologies, Inc. 2019 Printed in the USA, May 10, 2019 5994-0848EN

