

Agilent 850-DS Dissolution Sampling Station (G7930A) and Automated UV Systems Cary 50/60 (G7926A/G7927A) Supplies

Agilent Technologies is committed to optimizing your laboratory's productivity, so we have produced this list of the most commonly ordered supplies and parts for the Agilent 850-DS Dissolution Sampling Station and Automated UV Systems Cary 50/60.



Automated Sampling Replacement Tubings

Description	Comments	Part Number
Tubing assembly, sample and return cannula, 6-position	For 1 L 708-DS and 850-DS dissolution sampling station	1005-1920
Tubing assembly, sample and return cannula, 8-position		1005-1921
Tubing assembly, sample and return cannula, 6-position	For 2 L 708-DS and 850-DS dissolution sampling station	1005-1922
Tubing assembly, sample and return cannula, 8-position		1005-1923

Resident Sampling Kits

Description	Comments	Part Number
Resident sampling kit, 6-position, 900 mL	For 708-DS dissolution apparatus	K1001-01193
Resident sampling kit, 8-position, 900 mL		K1001-01194
Resident sampling kit, 6-position, 500 mL		K1001-01195
Resident sampling kit, 8-position, 500 mL		K1001-01196

UV Accessories

Description	Comments	Part Number
Flow cell only, 10 mm,	For Cary 50/60	11-1300
Flow cell only, 0.2 mm		11-1301
Flow cell only, 0.5 mm		11-1302
Flow cell only, 5 mm		11-1305
Flow cell only, 2 mm		11-1310
Flow cell only, 1 mm		11-1315
Tubing kit, dual path length	For Cary 60 fiber optic UV dissolution system	11-1320
Flow cell tubing replacement kit		11-1226
Fiber optic probe, 2 m	Without tip for 708-DS dissolution apparatus	11-1425
Fiber optic, 1 mm removable tip		11-1429
Fiber optic, 2 mm removable tip		11-1430
Fiber optic, 5 mm removable tip		11-1435
Fiber optic, 10 mm removable tip		11-1440
Fiber optic, 20 mm removable tip		11-1445
Fiber optic, 4 mm removable tip		11-1450

850-DS Sampling Vials, Tubes & Wellplates

Description	Comment	Part Number
A-Line screw top vial, 2 mL, clear, with write-on spot	100/pk	5190-9589
A-Line screw top vial, 2 mL, amber, with write-on spot	100/pk	5190-9590
A-Line crimp top vial, 2 mL, clear, with write-on spot	100/pk	5190-9591
A-Line crimp top vial, 2 mL, amber, with write-on spot	100/pk	5190-9592
Screw cap, bonded, blue, PTFE/white silicone septa	100/pk	5190-7021
Crimp cap, silver aluminum, PTFE/white silicone septa	100/pk	5182-0552
Glass tubes, 16x100 mm, 13 mL	250/pk	5022-6532
96-well plate, square wells, U shape, polypropylene, 2.2 mL	30/pk	5043-9300
96-sealing mat, square wells, pre-slit, silicone	50/pk	5043-9319

Thermal Printer Paper

Description	Comment	Part Number
Thermal printer paper	For use with 850-DS, 10 rolls/pk	G7910-87002

www.agilent.com/chem/dissolution

This information is subject to change without notice.