

# Agilent 400-DS Dissolution Apparatus 7 (G7975A) Supplies

## Apparatus 7—Small Volume Reciprocating Holder

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 400-DS Dissolution Apparatus.



### Dissolution Cells

Description	Part Number
Dissolution cell, 10 mL	33-9000
Dissolution cell, 5 mL	33-9005

### 400-DS Holder

Description	Part Number
PEEK basket	33-9042
PEEK basket assembly	33-9022
50-mesh basket, 10 mL	33-9029
50-mesh basket assembly, 10 mL	33-9041
Basket, mesh end, 20 m	33-9046
50-mesh basket, 5 mL	33-9051
50-mesh basket assembly, 5 mL	33-9052
Punctal plug holder, 5 mL	33-9047
Contact lens holder, 5 mL	33-9048
Contact lens holder, 16 mm, 5 mL	33-9050
Contact lens holder, 16 mm, 10 mL	33-9053
Flat blank holder, 5 mL	33-9049
Pacemaker lead holder, 1.7 mm, requires stent holder assembly	33-9025
Stent holder assembly, 5 mL	33-9009
Stent holder assembly, 10 mL	33-9012

### Sampling Vials

Description	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

[www.agilent.com/chem/dissolution](http://www.agilent.com/chem/dissolution)

This information is subject to change without notice.