

Supplies for Varian 705-DS Dissolution Apparatus

Agilent is committed to optimizing your laboratory's productivity. This reference guide shows the most commonly ordered supplies and parts for the Varian 705-DS Dissolution Apparatus. Keep this list so you can quickly find the supplies you need and minimize instrument downtime.



Description	Part Number
Basket Shafts/Assemblies	
Basket shaft, USP 3-clip, 21 inch	14-3620
Basket shaft, O-ring, 21 inch	14-3621
Basket shaft, PTFE-coated, 21 inch	14-3622
Basket shaft, USP 3-clip, 24 inch	14-3629
Basket shaft, for bolus basket, 24 inch	14-3630
Paddles	
Paddle, PTFE-coated, 21 inch	14-3594
Paddle, PTFE-coated, 24 inch	14-3596
Paddle, electropolished, 21 inch	14-3595
Paddle, electropolished, 24 inch	14-3597
Paddle, PEEK, 24 inch	14-3598
Mini paddle, PTFE-coated, 24 inch	14-3599
Mini paddle, electropolished, 24 inch	14-3600
Mega paddle, electropolished stainless steel, 24 inch	14-3601
Capsule Wires, Weights and Sinkers Baskets	
Capsule weight wire, 316 stainless steel, 0.032 in. dia., 50 ft.	12-3000
Capsule weight, 3-prong	12-3050
Capsule weight, robotic type, with magnetic cap	12-3055
Capsule weight, robotic type, with magnetic inserts in legs	12-3056
Sinker basket, Japanese Pharmacopoeia (JP)	12-3070

Description	Part Number
Baskets for Apparatus 1	
Basket, USP, 40-mesh	12-2100
Basket, USP, 40-mesh includes individual Certificate of Conformance	12-2100V
Basket, mini, 40-mesh	12-2102
Basket, gold-plated, 40-mesh	12-2105
Basket, PTFE-coated, 40-mesh	12-2110
Basket, 3-fin assembly, 40-mesh	12-2115
Basket, 10-mesh	12-2125
Basket, slotted, for suppository testing	12-2130
Basket, 325x2300-mesh, 5 µm	12-2146
Basket, 270x270-mesh, 53 µm	12-2151
Basket, 150x150-mesh, 104 µm	12-2152
Basket, 100x100-mesh, 140 µm	12-2154
Basket, 50x50-mesh, 280 µm	12-2155
Bolus basket, 20x20-mesh, 864 µm	12-2180
Bolus basket, 40x40-mesh, 381 µm	12-2185
Basket for Metformin, vertical sample holder, large	12-2165
Basket for Metformin, vertical sample holder, small	12-2170

Description	Part Number
Filters & Filter Tips for Sampling Cannulas	
Full Flow Filter, UHMWPE, blue, 10 µm, 1000/pk	17-4005
Full Flow Filter, UHMWPE, blue, 10 µm, 100/pk	17-4000
Full Flow Filter, UHMWPE, 1 µm, 100/pk	17-4003
Full Flow Filter, UHMWPE, 4 µm, 100/pk	17-4004
Full Flow Filter, UHMWPE, white, 35 µm, 1000/pk	17-4015
Full Flow Filter, UHMWPE, white, 35 µm, 100/pk	17-4010
Full Flow Filter, UHMWPE, red, 70 µm, 1000/pk	17-4025
Full Flow Filter, UHMWPE, red, 70 µm, 100/pk	17-4020
Full Flow Filter, PVDF, green, 10 µm, 1000/pk	17-4045
Full Flow Filter, PVDF, green, 10 µm, 100/pk	17-4040
Full Flow Filter, PVDF, yellow, 35 µm, 1000/pk	17-4055
Full Flow Filter, PVDF, yellow, 35 µm, 100/pk	17-4050
EaseAlign Vessels	
EaseAlign vessel, 1L	12-5000
EaseAlign vessel, 1L, with certificate	12-5000V
EaseAlign vessel, low actinic-red, 1L	12-5010
EaseAlign vessel, low actinic-red, 1L, with certificate	12-5010V
EaseAlign vessel, 100 mL	12-5040
EaseAlign vessel, 200 mL	12-5050
EaseAlign vessel, flat bottom for Enhancer Cell, 200 mL	12-5055
EaseAlign vessel, 2L	12-5070
EaseAlign vessel, 4L	12-5080
EaseAlign vessel, clear plastic, 1L	12-5200
EaseAlign Peak vessel, 1L	12-5500
EaseAlign Peak vessel, low actinic-red, 1L	12-5505
EaseAlign VF molded vessel, 1L	12-1502
EaseAlign VF molded vessel, 1L, with certificate	13-0020
EaseAlign Peak vessel, low actinic-red, 1L	12-5505
Vessel Conversion Kits and Accessories	
1L Conversion Kits	
Conversion kit, TruCenter, with ground groove, 1L	12-6100
Conversion kit, TruCenter, with manual sampling evaporation cover, 1L	12-6101
100 mL Conversion Kits	
Conversion kit, EaseAlign, 100 mL, without paddle	12-0310
200 mL Conversion Kits	
Conversion kit, EaseAlign, with 24 inch paddle, flat bottom vessel, 200 mL (for use with Enhancer Cells)	12-0301
Conversion kit, EaseAlign, 24 inch paddle, 200 mL vessel	12-0318
1L Conversion Kits	
Evaporation cover, EaseAlign, for use with 100/200 mL vessel	12-6315

Description	Part Number
Manual Sampling Accessories	
Bent cannula with adjustable gauge, 4.75 inch (900 mL)	12-3215
Bent cannula with adjustable gauge, 7.75 inch (900 mL)	12-3216
Adjustable gauge kit (cannulas not included)	12-3217
Sampling Cannulas and Accessories	
Resident sampling kit, 6-position, 900 mL	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
Centering Rings and Accessories	
Centering ring, EaseAlign, for 705-DS	12-6050
Locating pins, EaseAlign, 12/pk	12-6065
Alignment plate, TruCenter	12-6115
Intrinsic Dissolution Apparatus	
Intrinsic Dissolution Apparatus, 0.5 cm ² exposed surface area, with punch, shaft and holder	12-4100
Verification Tools	
Vessel centering tool, pass/fail gauge, EaseAlign centering rings, with certificate	12-7310V
Printer Paper	
Printer paper, 10 rolls/pk	12-9995
Printer paper, 100 rolls/pk	12-9996
Ribbon cartridge, 5/pk	12-9997

www.agilent.com/en/products/dissolution/apparatus

This information is subject to change without notice.

© Agilent Technologies, Inc. 2019
 Printed in the USA, June 28, 2019
 5994-1120EN