

Agilent 708-DS Dissolution Apparatus (6-position/G7910A and 8-position/G7911A) Supplies

Apparatus 1 – Rotating Basket Apparatus 2 – Rotating Paddles

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 708-DS Dissolution Apparatus.



Two-Piece Shafts for Apparatus 1

Description	Comment	Part Number
Upper receptor shaft, 21 in. (total length when assembled)	Two-piece basket shafts with individual certificate of conformance	14-3613
Receptor, 21 in. total, 6/pk	Convenient 6-pack	16-3613
Basket shaft, lower interchangeable, 3-clip, USP	Lower interchangeable basket shafts	14-3631
Basket, lower, 3-clip, 6/pk	Convenient 6-pack	16-3631
Basket shaft, lower interchangeable, O-ring	Lower interchangeable basket shafts	13-3632
Basket shaft, lower interchangeable, 3-clip, for 2 L		14-3634
Mini basket shaft, lower interchangeable, stainless steel, O-ring		13-3636

One-Piece Shafts for Apparatus 1

Description	Comment	Part Number
Basket shaft, USP 3-clip, 21 in.	With individual certificate of conformance	14-3620
Basket shaft, O-ring, 21 in.		13-3621
Basket shaft, PTFE-coated, 21 in.		13-3622
Basket shaft, USP 3-clip, 24 in.		13-3629
Basket shaft, for bolus basket, 24 in.		13-3630

Two-Piece Shafts for Apparatus 2

Description	Comment	Part Number
Upper receptor shaft, 21 in. (total length when assembled)	Two-piece paddle shaft with individual certificate of conformance	14-3613
Upper receptor shaft, for 15 in. (total length when assembled)		13-3612
Upper receptor shaft, 24 in. (total length when assembled)		13-3614
Receptor, 21 in. total, 6/pk	Convenient 6-pack	16-3613
Paddle, lower, electropolished stainless steel, 6/pk		16-3603
Paddle, lower, PTFE-coated, 6/pk	Convenient 6-pack	16-3602
Paddle	Lower interchangeable, PTFE-coated	14-3602
	Lower interchangeable, electropolished stainless steel 1 L	14-3603
	Lower interchangeable, PTFE-coated, for 2 L	13-3604
	Lower interchangeable, stainless steel, for 2 L	14-3605
	Lower interchangeable, PEEK	13-3606
Mini paddle	Lower interchangeable, PTFE-coated	14-3607
	Lower interchangeable, stainless steel	14-3608

One-Piece Paddle Shafts for Apparatus 2

Description	Part Number
Paddle, PTFE-coated, 21 in.	14-3594
Paddle, PTFE-coated, 24 in.	13-3596
Paddle, electropolished, 21 in.	14-3595
Paddle, electropolished, 24 in.	13-3597
Paddle, PEEK, 24 in.	13-3598
Mini paddle, PTFE-coated, 24 in.	14-3599
Mini paddle, electropolished, 24 in.	13-3600
Mega paddle, electropolished stainless steel, 24 in.	13-3601

Capsule Wires, Weights and Sinkers Baskets

Description	Part Number
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

Baskets for Apparatus 1

Description	Comment	Part Number
Basket, USP, 40-mesh		12-2100
Basket, USP, 40-mesh	With 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
	Vertical sample holder, small	12-2170

Shaft Accessories

Description	Comment	Part Number
Storage rack	For one-piece paddle/basket shafts	12-1350
Storage rack	for two-piece paddle/basket shafts	12-1352

Stationary Shafts and Basket

Description	Comment	Part Number
Stationary basket	Felodipine quadrangular	12-2069

Float-A-Lyzer Holder

Description	Comment	Part Number
Float-A-Lyzer holder	For use with USP Apparatus 1	12-2066
Float-A-Lyzer holder, stationary	For use with USP Apparatus 2	12-2067

Thermal Printer Paper

Description	Comment	Part Number
Thermal printer paper	For use with 708-DS, 10 rolls/pk	G7910-87002

Vessels and Conversion Kits

Description	Comment	Part Number
TruAlign VF molded vessel, clear, 1 L		13-0010
TruAlign vessel, 1 L, clear	With certificate	12-5149V
TruAlign vessel, low actinic - red, 1 L		12-5152V
TruAlign peak vessel, 1 L, clear		12-5153V
TruAlign peak vessel, low actinic - red, 1 L		12-5154
TruAlign vessel, 2 L, clear	With certificate	12-5157V
TruAlign vessel, low actinic - red, 2 L		12-5158V
TruAlign vessel, 100 mL		12-5145
TruAlign vessel, low actinic - red, 100 mL		12-5146
TruAlign vessel, 200 mL		12-5147
TruAlign vessel, low actinic - red, 200 mL		12-5148
TruAlign vessel, flat bottom, 200 mL (one per position)		12-5170
TruAlign vessel 100/200 mL conversion kit	Includes vessel, adapter and evaporation cover. Set of 6	12-6368
TruAlign vessel, 250 mL conversion kit, for Chinese Pharmacopoeia	Includes vessel, adapter, two-piece mini paddle, evaporation cover and manifold adapter. Set of 6	12-1450
TruAlign vessel, 250 mL conversion kit, for Chinese Pharmacopoeia	Includes vessel, adapter, two-piece mini paddle, evaporation cover and manifold adapter. Set of 8	12-1451

Evaporation Covers

Description	Comment	Part Number
708-DS evaporation cover, standard	for CP/USP sampling	K1005-05225
708-DS evaporation cover	low loss for use with resident sampling probes	K1005-05218

Manual Sampling Accessories

Description	Comment	Part Number
Bent cannula with luer lock	For sampling from 900 mL in 1 L vessels, 4.75 in.	12-3200
	For sampling from 500 mL in 1 L vessels, 7.75 in.	12-3210
Bent cannula with adjustable gauge	4.75 in. (900 mL), 708-DS	12-3221
	7.75 in. (500 mL), 708-DS	12-3222
Adjustable gauge kit, 708-DS		12-3223
Cannula	PEEK, for low-volume sampling	12-3219
Glass syringe, 20 cc		12-3230

Full Flow Filters

Description	Comment	Part Number	
Full flow filter	UHMWPE, blue, 10 µm, 1000/pk	17-4005	
	UHMWPE, blue, 10 µm, 100/pk	17-4000	
	UHMWPE, 1 µm, 100/pk	17-4003	
	UHMWPE, 4 µm, 100/pk	17-4004	
	UHMWPE, white, 35 µm, 1000/pk	17-4015	
	UHMWPE, white, 35 µm, 100/pk	17-4010	
	UHMWPE, red, 70 µm, 1000/pk	17-4025	
	UHMWPE, red, 70 µm, 100/pk	17-4020	
	PVDF, green, 10 µm, 1000/pk	17-4045	
	PVDF, green, 10 µm, 100/pk	17-4040	
	PVDF, yellow, 35 µm, 1000/pk	17-4055	
	PVDF, yellow, 35 µm, 100/pk	17-4050	
	Filter tip	PE, 2 µm, 100/pk	17-4100
		PE, 5 µm, 100/pk	17-4110
PE, 10 µm, 100/pk		17-4120	

www.agilent.com/chem/dissolution

This information is subject to change without notice.

© Agilent Technologies, Inc. 2018
Printed in the USA, November 1, 2018
5994-0388EN