

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

## Apparatus 5—Paddle over Disk Apparatus 6—Rotating Cylinder

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.



### Upper Shafts for Apparatus 5 and 6

| Description   | Comment   | Part Number |
|---|---|-------------|
| Upper receptor shaft, 21 in.<br>(total length when assembled)     | Two-piece paddle shaft with individual certificate of conformance | 14-3613     |
| Upper receptor shaft, for 15 in.<br>(total length when assembled) |   | 13-3612     |
| Upper receptor shaft, 24 in.<br>(total length when assembled)     |   | 13-3614     |

### Lower Interchangeable Two-Piece Shafts for Apparatus 5

| Description | Comment   | Part Number |
|-------------|---|-------------|
| Paddle      | Lower interchangeable, PTFE-coated                        | 14-3602     |
| Paddle      | Lower interchangeable, electropolished stainless steel 1L | 14-3603     |
| Paddle      | Lower interchangeable, PEEK                               | 13-3606     |
| Mini paddle | Lower interchangeable, PTFE-coated                        | 13-3607     |
| Mini paddle | Lower interchangeable, stainless steel                    | 14-3608     |

### One-Piece Paddle Shafts for Apparatus 5

| Description                          | Comment                                    | Part Number |
|--------------------------------------|--|-------------|
| Paddle, PTFE-coated, 21 in.          | With individual certificate of conformance | 13-3594     |
| 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     |

## Paddle over Disk

| Description  | Part Number |
|--|-------------|
| Paddle over disk assembly, with 35 mm opening, 40-mesh screen  | 12-4200     |
| Paddle over disk assembly, 125 µm, EP                          | 12-4201     |
| Replacement screens, 125 µm, EP, 100/pk                        | 12-4209     |
| Replacement screens, 35 mm openings, 40-mesh, 100/pk           | 12-4210     |
| Disk disassembly kit, for 12-4200                              | 12-4211     |
| Paddle over disk assembly, with 56 mm opening, 120-mesh screen | 12-4230     |
| Replacement screens, 56 mm openings, 120-mesh, 100/pk          | 12-4231     |
| Disk disassembly kit, for 12-4230                              | 12-4232     |

## Transdermal Patch Holders and Accessories

| Description                                    | Part Number |
|--|-------------|
| Transdermal patch holder, 2.5 cm <sup>2</sup>  | 12-4300     |
| Transdermal patch holder, 5.0 cm <sup>2</sup>  | 12-4310     |
| Transdermal patch holder, 7.0 cm <sup>2</sup>  | 12-4320     |
| Transdermal patch holder, 10.0 cm <sup>2</sup> | 12-4330     |
| Cuprophane membrane, 126x345 mm, 10 sheets/pk  | 12-1370     |

## Suspension Cups

| Description  | Part Number |
|--|-------------|
| Weighted holder, 5.72 cm outside dia., 0.52 cm inside dia. | 12-4050     |
| Insert cup, 17.5x8.2 mm, use with 12-4050                  | 12-4055     |
| Insert cup, 20x8.2 mm, use with 12-4050                    | 12-4060     |

## Rotating Cylinder and Accessories

| Description  | Part Number |
|--|-------------|
| Lower interchangeable shaft for rotating cylinder                                    | 12-1371     |
| Lower interchangeable shaft for rotating cylinder (with certificate)                 | 12-1371V    |
| Rotating cylinder, 25 mm outside dia., with interchangeable shaft                    | 12-1372     |
| Rotating cylinder, 25 mm outside dia., with interchangeable shaft (with certificate) | 12-1372V    |
| Height gauge, rotating cylinder, USP   | 12-7335     |
| Cuprophane membrane, 126x345 mm, 10 sheets/pk  | 12-1370     |

## Intrinsic Dissolution and Accessories

| Description   | Part Number |
|---|-------------|
| Intrinsic dissolution apparatus, 0.5 cm <sup>2</sup> exposed surface area, with punch, shaft and holder, for 7000E/7010, 708-DS, 705-DS | 12-4101     |
| Intrinsic dissolution apparatus, 0.125 cm <sup>2</sup> exposed surface area (18 in. shaft)  | 12-4110     |
| Intrinsic die, 0.5 cm <sup>2</sup> exposed surface area   | 12-4120     |
| Punch   | 12-4140     |
| Shaft and die holder only, for intrinsic dissolution (18 in. shaft)   | 12-4150     |
| Surface plate, for intrinsic dissolution  | 12-4130     |

## 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     |

## Vessels

| Description                                  | Comment          | Part Number |
|--|------------------|-------------|
| TruAlign VF molded vessel, clear, 1 L        | With certificate | 13-0010     |
| TruAlign vessel, 1 L, clear                  | With certificate | 12-5149V    |
| TruAlign vessel, low actinic - red, 1 L      | With certificate | 12-5152V    |
| TruAlign peak vessel, 1 L, clear             | With certificate | 12-5153V    |
| TruAlign peak vessel, low actinic - red, 1 L |                  | 12-5154     |

## Enhancer Cell

| Description   | Part Number |
|---|-------------|
| Enhancer cell, 4 cm <sup>2</sup> surface area membrane (one per position)                 | 12-4000     |
| Enhancer cell, 2 cm <sup>2</sup> surface area membrane (one per position)                 | 12-4001     |
| Enhancer cell, 0.5 cm <sup>2</sup> surface area membrane (one per position)               | 12-4002     |
| Height spacer (1 cm)/alignment tool, for 4 cm <sup>2</sup> enhancer cell (one required)   | 12-4020     |
| Height spacer (1 cm)/alignment tool, for 2 cm <sup>2</sup> enhancer cell (one required)   | 12-4021     |
| Height spacer (1 cm)/alignment tool, for 0.5 cm <sup>2</sup> enhancer cell (one required) | 12-4022     |
| Adjustment tool, for all enhancer cell sizes (one required)                               | 12-4015     |

## 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     |
| Bent cannula with luer lock        | 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     |
| Bent cannula with adjustable gauge | 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     |

## Filters and Filter Tips for Sampling Cannulas

| Description      | Comment                       | Part Number |
|------------------|-------------------------------|-------------|
| 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     |
| Filter tip       | PE, 2 µm, 100/pk              | 17-4100     |
| Filter tip       | PE, 5 µm, 100/pk              | 17-4110     |
| Filter tip       | PE, 10 µm, 100/pk             | 17-4120     |

## Thermal Printer Paper

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

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

This information is subject to change without notice.

© Agilent Technologies, Inc. 2018  
Printed in the USA, November 1, 2018  
5994-0392EN