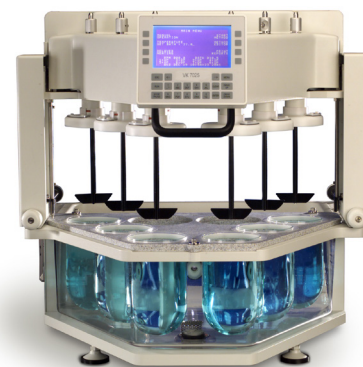


## Supplies for Varian VK7020/7025 and 7030 V-Series Dissolution Testers

Agilent is committed to optimizing your laboratory's productivity. This reference guide shows the most commonly ordered supplies and parts for the Varian VK7020/7025 and 7030 V-Series Dissolution Testers. Keep this list so you can quickly find the supplies you need and minimize instrument downtime.



Description	Part Number
<b>Basket Shafts/Assemblies</b>	
Basket shaft 14.5 inch (37 cm), USP	14-3615
<b>Paddles</b>	
Paddle, PTFE-coated 14.5 inch (37 cm)	14-3589
Paddle, electropolished 14.5 inch (37 cm)	14-3590
Paddle, PEEK 14.5 inch (37 cm)	14-3591
Paddle, electropolished, 200 mL 14.5 inch (37 cm)	14-3593
<b>Capsule Wires, Weights and Sinkers Baskets</b>	
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
<b>Baskets for Apparatus 1</b>	
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
<b>Filters &amp; Filter Tips for Sampling Cannulas</b>	
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
<b>Replacement Sampling Accessories</b>	
Sample tubing/cannulas, sampling/return, 8-position	1005-1786
Sample tubing/cannulas, sampling/return, 6-position	1005-1787
<b>TruCenter Vessels</b>	
TruCenter vessel, without adapter, 200 mL	12-5033
TruCenter vessel, with adapter, 200 mL	12-5034
TruCenter vessel, with collar, 1L	12-5035
TruCenter vessel, with collar, 1L, with certificate	12-5035V
TruCenter vessel, low actinic-red, with collar, 1L	12-5120
TruCenter vessel, low actinic-red, with collar, 1L, with certificate	12-5120V
TruCenter Peak vessel, with collar, 1L	12-5125
TruCenter VF molded vessel, 1L	12-1503
TruCenter VF molded vessel, 1L, with certificate	13-0030
<b>Conversion Kits and Accessories</b>	
<b>200 mL Conversion Kit</b>	
TruCenter conversion kit, with paddle, 200 mL	12-0312
TruCenter conversion kit, with paddle and temperature monitoring, 200 mL	12-0314
<b>Conversion Kits Accessories</b>	
TruCenter vessel, with adapter, 200 mL	12-5034
TruCenter vessel, without adapter, 200 mL	12-5033

Description	Part Number
<b>Evaporation Covers and Accessories</b>	
Evaporation cover, manual sampling, opening for Japanese sinkers	12-6352
Evaporation cover, dual sampling	12-6353
<b>Printer Paper</b>	
Printer paper, 10 rolls/pk	12-9995
Printer paper, 100 rolls/pk	12-9996
Ribbon cartridge, 5/pk	12-9997
<b>Intrinsic Dissolution Apparatus</b>	
Intrinsic Dissolution Apparatus, 0.5 cm <sup>2</sup> exposed surface area	12-4105
Intrinsic Dissolution Apparatus, 0.125 cm <sup>2</sup> exposed surface area	12-4111

[www.agilent.com/en/products/dissolution/apparatus](http://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  
 5994-1122EN