

Level 1 of VANQUISH Sample Container Kit

Kit catalogue number: **VSCK**

Vanquish Sample Container Kit

The Vanquish system offers a wide variety of supported vials and well plates. This sample container gives you an overview of the recommended and proven vial options in line with the required sample racks.

Microsampling Vials (6 mm OD):

They have a usable sample volume of 2 - 250 µL. The dedicated sample rack for the 6 mm OD vials provides 96 positions. This rack is optional.

Standard & Reduced Volume Vials (12 mm OD):

Available as screw, crimp or snap cap closure systems. They have a usable volume of 2 - 1500 µL depending on the vial type aligned to your sample amount requirements. Single micro inserts can be used for volume reduction down to 2 - 200 µL as well. The dedicated sample rack for the 12 mm OD vials has 54 positions. This sample rack is shipped with the Vanquish autosampler.

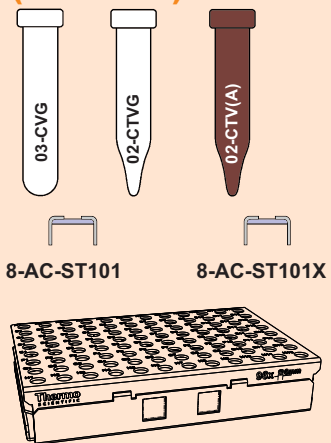
Increased Volume Vials (15 mm OD):

These vials have a usable volume of 0.5 - 3.5 mL. The sample rack can be ordered as accessory. (Please contact your local Thermo Fisher Scientific sales organization for HPLC products)

Large Volume Vials (22.5 mm OD):

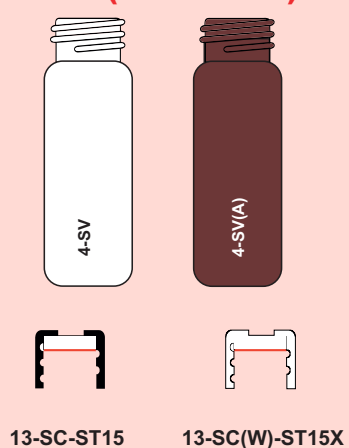
The Vanquish autosampler has the capacity of up to 12 x 22.5 mm OD vials in its carousel.

Microsampling Vials (6 mm OD)

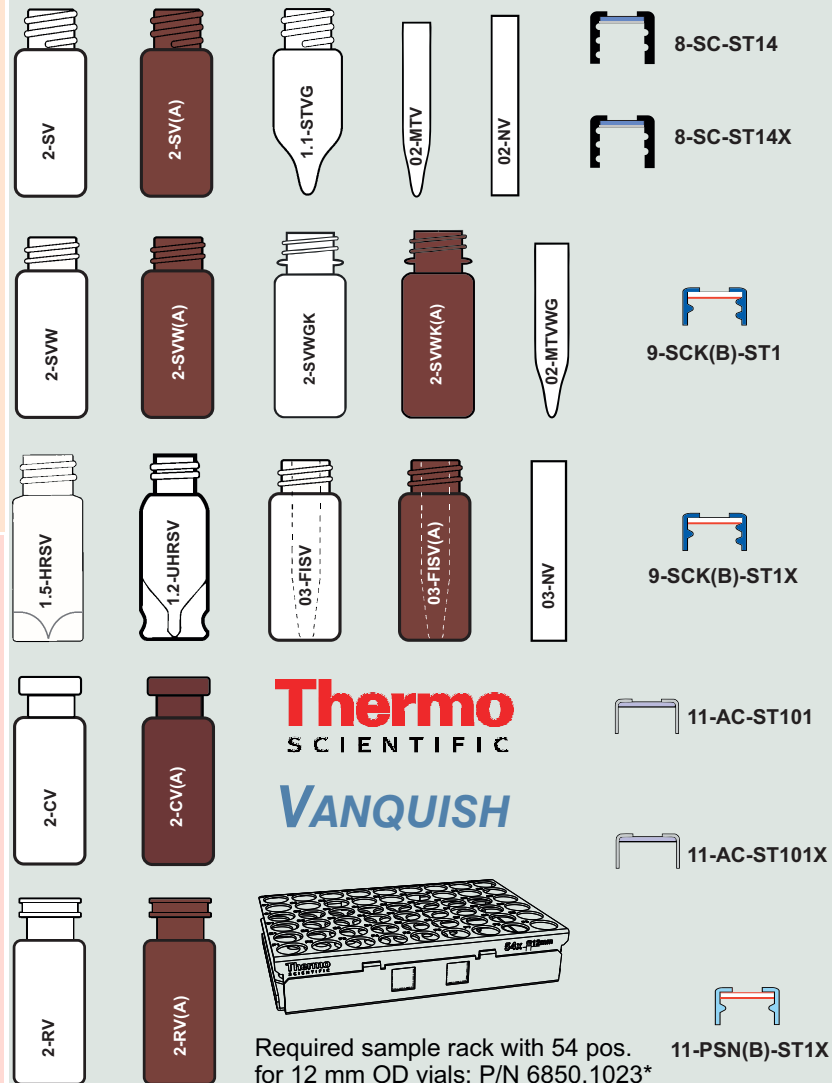


Required sample rack with 96 pos. for 6 mm OD vials: P/N: 6850.1026*

Increased Volume Vials (15 mm OD)

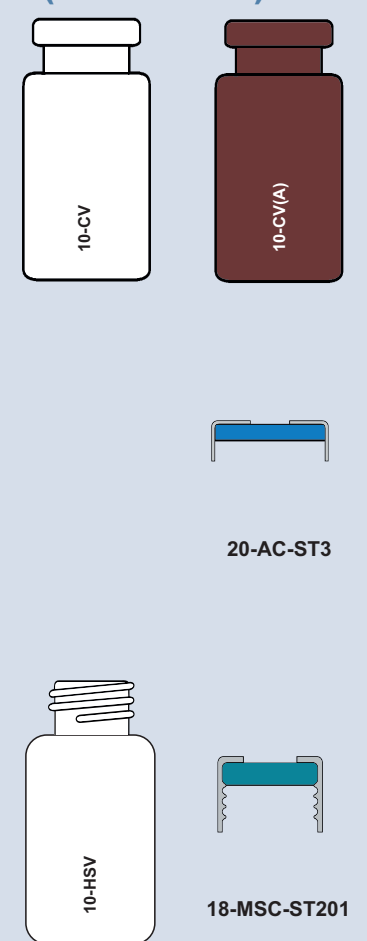


Standard & Reduced Volume Vials (12 mm OD) & Inserts



Required sample rack with 54 pos. for 12 mm OD vials: P/N 6850.1023*

Large Volume Vials (22.5 mm OD)



© 2014 Thermo Fisher Scientific Inc. All rights reserved.
All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change.
Not all products are available in all countries. Please consult your local sales representative for details.

For more vials& closures information visit www.thermoscientific.com/chromatography

*For more information about instrument accessories visit www.thermoscientific.com

USA and Canada +1 800 332 3331
France +33 (0)1 60 92 48 34
Germany +49 (0) 2423 9431 20 or 21
United Kingdom +44 (0)1928 534110
Japan +81 3 5826 1615

China +86 21 68654588 +86 10 84193588
+86 20 83145199 800 810 5118
India +91 22 6742 9494 +91 27 1766 2352
Australia 1 300 735 292 (free call domestic)
New Zealand 0800 933 966 (free call domestic)
All Other Enquiries +44 (0) 1928 534 050

Technical Support
North America +1 800 332 3331
Outside North America +44 (0) 1928 534 440

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand

Level 2 of VANQUISH Sample Container Kit

Kit catalogue number: **VSCK**

Vanquish Sample Container Kit

The Vanquish system offers you a huge variety of supported vials and well plates. This content sheet gives you an overview of the recommended and proven well plate & seal options.

No adapter or rack is required for the usage of well plates in the Vanquish system.

384 Square Deep Well Plate for volumes from 5 - 240 μ L, Chromatography CERTIFIED

96 Round Deep Well Plate for volumes from 50 - 900 μ L, Chromatography CERTIFIED

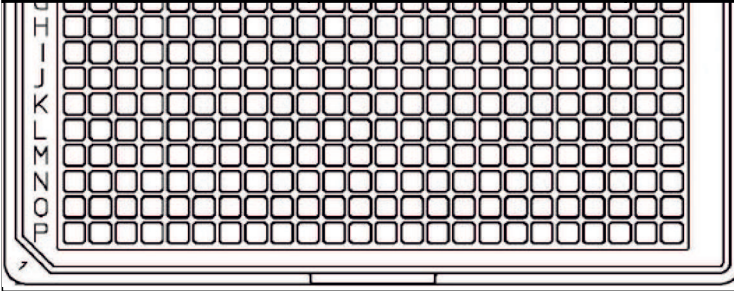
96 Round Shallow Well Plate for volumes from 10 - 190 μ L, Chromatography CERTIFIED

384 Square Well Silicone Mat, Pre-slit

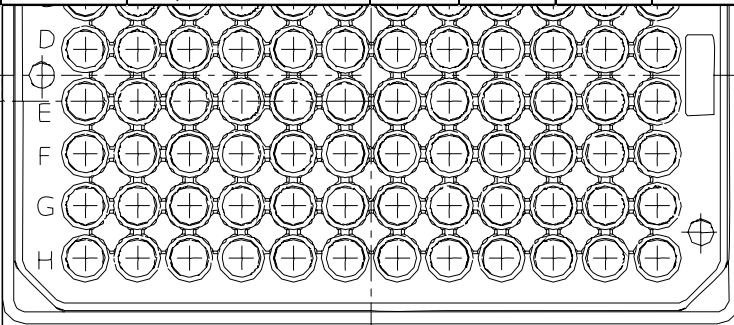
96 Round Well Silicone Mat, Pre-slit

96 pre-slitted and resealing Tape with no glue at the cavity area, polar organic solvent stable.

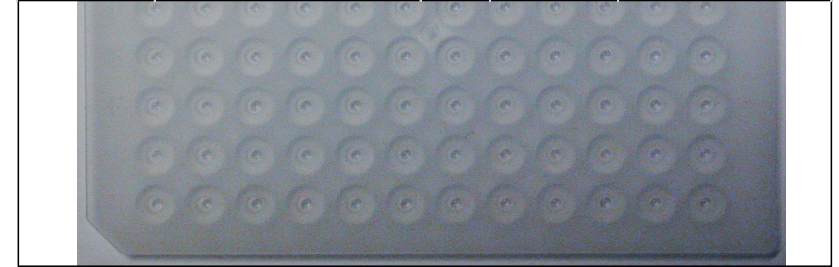
Cat. No.	Description	Total Height (mm)	Well Format	Total Volume (μ L)	working volume range μ L/well
60180-P204	PP 384 Square Deep Well Plate, CERTIFIED	30.2	V-Shape, square	300	5 - 240



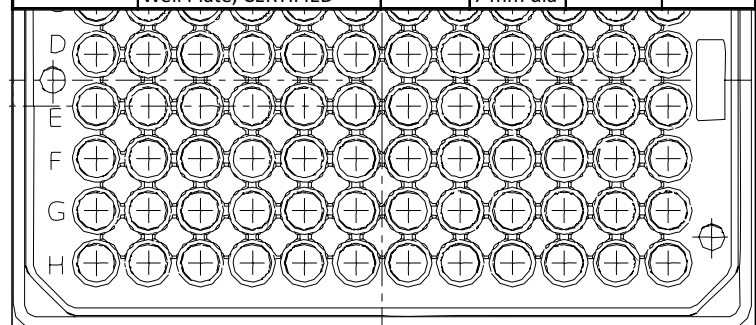
Cat. No.	Description	Total Height (mm)	Well Format	Total Volume (μ L)	working volume range μ L/well
60180-P201	PP 96 Round Deep Well Plate, CERTIFIED	41.6	U-Shape, 7 mm dia	1000	50 - 900



Cat. No.	Description	Color	Material	Well Design
60180-M113	MicroMat CLR Silicone Mat	Clear	Silicone	96 Round Well – Pre-slit, 7 mm dia
60180-M150	MicroMat CLR Silicone Mat	Clear	Silicone	384 Square Well – Pre-slit
60180-M146	3 layer sealing tape, no adhesive at cavity area	Clear	PET/Silicone/PET	Pre-slit



Cat. No.	Description	Total Height (mm)	Well Format	Total Volume (μ L)	working volume range μ L/well
60180-P207	PP 96 Round Shallow Well Plate, CERTIFIED	14.7	V-Shape, 7 mm dia	220	10 - 190



© 2014 Thermo Fisher Scientific Inc. All rights reserved.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

For more information about well plates visit www.thermoscientific.com/webseal

USA and Canada +1 800 332 3331
France +33 (0)1 60 92 48 34
Germany +49 (0) 2423 9431 20 or 21
United Kingdom +44 (0)1928 534110
Japan +81 3 5826 1615

China +86 21 68654588 +86 10 84193588
 +86 20 83145199 800 810 5118
India +91 22 6742 9494 +91 27 1766 2352
Australia 1 300 735 292 (free call domestic)
New Zealand 0800 933 966 (free call domestic)
All Other Enquiries +44 (0) 1928 534 050

Technical Support
North America +1 800 332 3331
Outside North America +44 (0) 1928 534 440

Thermo
 SCIENTIFIC

A Thermo Fisher Scientific Brand