

# GoPure™ 96-Well Screening Plates

Pub. No. MAN1000117 Rev. A00



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

## Product description

POROS™ GoPure™ 96-Well Screening Plates and CaptureSelect™ GoPure™ 96-Well Screening Plates allow for fast screening of multiple chromatographic conditions in parallel.

**Table 1 Resin characteristics**

Resin	Chemistry	Surface functionality
AAVX	Affinity	AAVX ligand
AAV9	Affinity	AAV9 ligand
Oglio (dT)25	Affinity	Poly-deoxythymidine (dT25-mer) ligand
MabCaptureC	Affinity	Protein A ligand (Manufactured by Thermo Fisher Scientific)
FcXP	Affinity	VHH affinity ligand

## Specifications

**Table 2 Well plate specifications**

Condition	Specification
Resin volume per well	20 µL
Buffer solution	18-20% (v/v) aqueous ethanol
Resin slurry volume per well	220 µL
Well volume	800 µL
Plate material	Polypropylene
Frit	25 µm polyethylene filter frit
Storage temperature	2-8°C Do not freeze.

## Ordering information

Product	Resin volume per well	Cat. No.
POROS™ GoPure™ AAVX 96-Well Screening Plates	20 µL	A65809
POROS™ GoPure™ AAV9 96-Well Screening Plates	20 µL	A65810
POROS™ GoPure™ Oligo (dT)25 96-Well Screening Plates	20 µL	A65811
GoPure™ MabCaptureC™ 96-Well Screening Plates	20 µL	A65812
CaptureSelect™ GoPure™ FcXP 96-Well Screening Plates	20 µL	A65813

## Related documentation

Publication	Pub. No.
<i>CaptureSelect™ GoPure™ 96-Well Screening Plates User Guide for plasma applications</i>	MAN1000118
<i>GoPure™ MabCaptureC™ 96-Well Screening Plates User Guide for protein antibody applications</i>	MAN1000119
<i>POROS™ GoPure™ 96-Well Screening Plates User Guide for gene therapy applications</i>	MAN1000120
<i>POROS™ GoPure™ 96-Well Screening Plates User Guide for mRNA purification applications</i>	MAN1000121

## Support

For service and technical support, go to [thermofisher.com/poros](https://www.thermofisher.com/poros) or call toll-free in US: 1.800.955.6288.

For the latest service and support information at all locations, or to obtain Certificates of Analysis or Safety Data Sheets (SDSs; also known as MSDSs), go to [thermofisher.com/support](https://www.thermofisher.com/support) or contact your local Thermo Fisher Scientific representative.

## Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](https://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](https://www.thermofisher.com/support).



**Revision history:** Pub. No. MAN1000117 A00

Revision	Date	Description
A00	1 May 2024	New document for the release of POROS™ GoPure™ AAVX 96-Well Screening Plates (Cat. No. A65809), POROS™ GoPure™ AAV9 96-Well Screening Plates (Pub. No. A65810), POROS™ GoPure™ Oligo (dT)25 96-Well Screening Plates (Cat. No. A65811), GoPure™ MabCaptureC™ 96-Well Screening Plates (Cat. No. A65812), and CaptureSelect™ GoPure™ FcXP 96-Well Screening Plates (Cat. No. A65813).

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.