# **INSTRUCTIONS**

# POROS XQ and HQ Resins

# 82073 82074 82077 82078

| Number | Description                     |
|--------|---------------------------------|
| 82073  | POROS XQ, 10mL settled resin    |
| 82074  | POROS XQ, 50mL settled resin    |
| 82077  | POROS 50 HQ, 10mL settled resin |
| 82078  | POROS 50 HQ, 50mL settled resin |

Resin: Crosslinked poly(styrene-divinylbenzene) Supplied: 50% slurry in 20% ethanol

Storage: Upon receipt store at 2-30°C. DO NOT FREEZE. Product shipped on wet ice.

# **Table of Contents**

| ntroduction                                   | 1 |
|---|---|
| mportant product information                  | 2 |
| Resin handling                                | 2 |
| Column packing for lab-scale columns (<100mL) | 3 |
| Qualify the column                            | 4 |
| Purification: Capture and elute               | 5 |
| Purification: Flow-through/Polishing          | 6 |
| Cleaning-in-place and resin storage           | 7 |
| Froubleshooting                               | 7 |
| Related Thermo Scientific Products            | 9 |
|   |   |

# Introduction

Thermo Scientific<sup>TM</sup> POROS<sup>TM</sup> Strong Anion Exchange Resins (POROS AEX resins) are designed for charge-based chromatographic separation of biomolecules including recombinant proteins, monoclonal antibodies, DNA, viruses, and peptides over a wide range of pH, conductivity and flow conditions. Two different surface chemistries are available with each chemistry offering unique selectivity that can impact both capacity and separation in a process.

Thermo Scientific<sup>TM</sup> POROS<sup>TM</sup> XQ resin is functionalized with fully quaternized quaternary amine. Thermo Scientific<sup>TM</sup> POROS<sup>TM</sup> 50 HQ resin is functionalized with quaternized polyethyleneimine groups (strong with some weak AEX functionality). The salt tolerance of POROS XQ resin typically allows proteins to be loaded and bound to the resin in buffer solutions that may limit binding to other supports, such as those containing  $\leq 150$ mM NaCl. POROS XQ resin provides a higher binding capacity with higher salt tolerance when compared to Thermo Scientific<sup>TM</sup> POROS<sup>TM</sup> 50 HQ resin.

The resin backbone of both POROS XQ and 50 HQ consists of a rigid, highly porous, crosslinked poly[styrene divinylbenzene] with an additional polyhydroxyl surface coating to ensure low nonspecific binding. The intrinsic stability and open pore structure of POROS beads allow rapid binding of large molecules, consistently producing high dynamic binding capacities over a wide range of flow conditions. POROS resins improve the speed, flexibility, and optimization of purifications when compared to soft supports (e.g., agarose), thus increasing performance of purifications over a wide range laboratory and process-scale applications.



| Table 1. Properties of Ther | mo Scientific POROS 50 HQ and XQ Resins.   |
|-----------------------------|--|
| Support matrix              | Crosslinked poly(styrene-divinylbenzene)   |
| Surface functionality       | 50 HQ: Quarternized polyethyleneimine (Strong)   |
|                             | XQ: Proprietary quaternary amine (Strong)  |
| Dynamic binding capacity    | 50 HQ: 76-106mg/mL <sup>[1]</sup>  |
|                             | XQ: ≥140mg/mL <sup>[2]</sup>   |
| Shipping solvent            | 20% ethanol  |
| Average particle size       | 50µm   |
| Shrinking/swelling          | <1% from 1-100% solvent  |
| Mechanical resistance       | 100bar/1450psi/10MPa   |
|                             | Compatibility  |
| pH range                    | 1-14   |
| Ionic strength range        | 0-5M, all common salts   |
| Buffer additives            | All common reagents including 1M sodium hydroxide, 8M urea,<br>6M guanidine hydrochloride, ethylene glycol, and detergents   |
| Solvents                    | Water, 0-100% alcohol, acetonitrile, 1-2M acids (e.g., acetic, hydrochloric, phosphoric), and other common organic solvents  |
|                             | Do not expose resin to strong oxidizers (e.g., hypochlorite), oxidizing acids (e.g., nitric acid), strong reducing agents (e.g., sulfite), acetate, or benzyl alcohol. |
| Operating temperature       | 2-30° C<br>Do not freeze   |

1] 50% breakthrough of BSA in 20mM tris, pH 8.0 at 100cm/hr in 0.46cm ID x 10cm length column

[2] 5% breakthrough of BSA in 20mM tris, 50mM NaCl, pH 8.0 at 300cm/hr in 0.46cmD x 20cm length column

#### Important product information

- In the presence of low-conductivity solutions, POROS XQ resins have electrostatic properties that may cause the resin to stick to glass and some types of plastic. To aid in resin handling, resin can be adjusted into a slurry containing 0.1M NaCl.
- For all column chromatography applications we recommend that highly pure buffer components be used and that all buffers and samples are degassed and filtered through a 0.45µm filter before use.
- Final protein yield and purity during a chromatographic separation are dependent upon the sample complexity and concentration in load sample, as well as the buffer conditions and flow rates used. Therefore, it is important to optimize these parameters before attempting a large-scale purification.
- If you are required to stir the resin slurry over an extended period of time, use a free-standing paddle stirrer. Do not use magnetic stir bars or paddles that mate with the bottom of the container.

# Resin handling (<100mL)

Below are general recommendations for handling resin, adjusting slurry percentage, and exchanging resin into new buffers.

- 1. Transfer the required volume of resin slurry to a bottle-top filter or sintered-glass frit.
- 2. Apply vacuum to remove the liquid from the top of the resin bed.
- 3. Resuspend the resin cake to the desired final resin slurry volume with 0.1M NaCl. Mix with a plastic or rubber spatula. Do not grind or press resin against the sintered glass material. If using a paper filter membrane, take care to not damage or rip the filter material.
- 4. Repeat steps 2 and 3 for a total of three exchanges. After the third exchange, recover the resin slurry into an appropriate container and store appropriately.
- 5. Verify the slurry concentration by transferring 10-100mL of slurry into a graduated cylinder and allow the resin to settle for  $\geq$ 4 hours. If needed, adjust the slurry concentration to 50–70%.
- 6. Proceed with column packing (below).



# Column packing for lab-scale columns (<100mL)

**Note:** POROS resins are mechanically rigid and incompressible and can be packed effectively in low-pressure glass columns and high-pressure stainless steel columns. We do not recommend packing POROS XQ in water, ethanol, or low-conductivity solutions because of the high ionic surface charge of POROS anion exchange resin. For volumes greater than 100mL, please refer to the POROS column packing instructions at https://tools.thermofisher.com/content/sfs/manuals/4468731.pdf.

There is a 1.08 packing factor for POROS XQ resin and a 1.06 packing factor for POROS HQ resins to account for the difference in bed volume between a gravity-settled bed in 0.1M NaCl and a 1- to 3-bar pressure-packed bed. This factor, along with the slurry ratio, is used to determine the volume of slurry required to yield the intended final column volume.

1. Determine the required slurry volume:

 $Required \ slurry \ volume = \frac{Target \ Column \ Volume}{Slurry \ ratio} \times packing \ factor$ 

Example: A column of dimensions 2.5cm ID × 10.2cm length would have a packed bed of 50mL (*Volume* =  $\pi r^2 H$ )

To properly pack this volume using a 50% resin slurry of POROS 50 HQ, 106mL is needed:

$$\frac{50mL}{0.50} \times 1.06 = 106mL$$
 of slurry required

- 2. Ensure that the column outlet is closed and plumbed directly to waste. Do not connect the column outlet to the chromatography system. Plumbing into the system creates backpressure that fights against the inlet pressure trying to settle the bed and pack the column.
- 3. Ensure that the column is level and locked in place before beginning the pack. If needed, attach appropriate column packing extender to allow full resin slurry volume to be added.
- 4. Add 0.1M NaCl to the empty column until the solution is 5cm above the bottom frit.
- 5. Wet bottom frit by opening the column outlet and allow liquid to enter frit and begin to flow. This will ensure the air trapped into the frit will not enter resin bed. Close outlet when the solution reaches 3-5mM above the bottom frit
- 6. Slowly deliver the required slurry volume to the column and immediately move on to the next steps. Do not allow the resin to fully gravity-settle in the column before packing.

Note: POROS resin beads typically have a density similar to water and do not settle rapidly.

- 7. Bring the top flow adaptor to the top of the slurry level and tighten the O-ring until it begins to form a seal with the column. Do not push the resin up and over the O-ring.
- 8. Continue to lower the adaptor slowly to remove the bubbles from the top of the slurry liquid. Do not allow large air bubbles between the top adaptor and the top of the resin slurry.
- 9. Tighten the O-ring completely and connect the column to the system.
- 10. Open the column outlet.
- 11. Start flow and increase the flow rate to  $\geq$ 50% above the maximum flow rate used during your planned purifications. For columns with a diameter of <1.0cm, a flow rate of 1000-2000cm/hr is recommended.
- 12. Continue flow until a stable resin bed forms. Monitor the pressure; it will gradually rise as the column packs. Do not exceed maximum pressure recommended for column or instrument hardware.

**Note:** While packing the column, you may observe some turbidity in the eluent as packing begins. Turbidity will clear as packing proceeds and 1-2 bed volumes of packing buffer pass through the column.

13. After the bed is formed, remove the top adaptor and remove excess solution with a pipette. Leave 2-5cm of liquid above the resin bed.

Thermofisher.com



- 14. Re-insert top adaptor and bring the top flow adaptor to the top of the resin bed. Tighten the O-ring until it begins to make a seal. Lower the adaptor slowly to remove the bubbles from the top of the column.
- 15. Bring the adaptor into contact with the top of the bed and seal O-ring. POROS resin does not shrink or swell, so an open headspace is not recommended. Do not push the resin over the O-ring or use pressure to compress the resin bed.
- 16. Flow at the packing flow rate again for 2-5 column volumes, taking note of the bed height and pressure at the desired flow rate.
- 17. If running the column in a unidirectional flow, packing is completed. If running the column in both directions, continue on to the next step to further condition the column.
- 18. If reversing the flow of the column during operation, condition the column in upflow. Switch column inlet and outlet position related to the pump flow for upflow and downflow. Do not invert the column.
  - a. Flow 2–3 column volumes in upflow at the operating flow rate.
  - b. Flow 2–3 column volumes in downflow at the operating flow rate, then adjust the adaptor if needed.
  - c. Flow 2 column volumes after you adjust the adaptor.

### Qualify the column

To qualify the integrity of a packed column, determine the HETP (height equivalent to a theoretical plate) and asymmetry using a non-binding analyte (a "plug"). Equation calculators are available in the Thermo Scientific<sup>™</sup> ChromaPro<sup>™</sup> mobile application available online. Alternatively, use the equations following the protocol.

#### A. Setting specifications

Qualification results depend on a number of factors including:

- Solutions and method used
- Scale
- Column hardware
- Chromatography system

After defining a column qualification procedure for a specific system (column plus chromatography system), base the qualification acceptance criteria on historical values and ranges instead of theoretical qualification results. Performing the column qualification method consistently and reproducibly is critical to obtaining meaningful results.

| Condition            | Recommendation                       |                                      |
|----------------------|--------------------------------------|--------------------------------------|
| Resin                | XQ                                   | 50 HQ                                |
| Flow rate            | Target operating flow rate (cm/hour) | Target operating flow rate (cm/hour) |
| Equilibration buffer | 0.5M NaCl                            | 0.1M NaCl                            |
| Plug solution        | 2M NaCl                              | 1M NaCl                              |
| Plug volume          | 4% of column volume                  | 2% of column volume                  |

#### Table 2. Recommended column qualification conditions.

#### **B.** Qualification protocol

- 1. Equilibrate the column and all buffers to working temperature. Ensure that all solutions are degassed.
- 2. Equilibrate the system.
- 3. Equilibrate the column with 5-15 column volumes of the equilibration buffer.
- 4. Inject your plug and continue running the system for 1.5 column volumes.
- 5. Record the elution volume peak shape of the plug after it passes through the column
- 6. Determine HETP and Asymmetry factor.







# **Purification: Bind and elute**

In a bind and elute method, sample components are bound to the resin and then eluted with increasing salt concentration or modification of pH conditions to sequentially elute sample components at different conditions. When performing IEX chromatography some optimization is always required to get optimal separation. The most common areas to optimize include the buffer conditions used for binding and elution, as well as the number of column volumes over which an elution gradient is applied. The following notes and protocol are provided as a starting point. Adjust as appropriate for your sample and the desired level of separation required. See the Troubleshooting section for more tips on optimizing protocols.

| Buffer composition  | • To ensure resin binding, the binding buffer/sample buffer pH should be 1 to 3 units above the pI of the target molecule. There is a direct correlation between dynamic binding capacity and the difference between the pI and pH of your buffer. Dynamic binding capacity will increase as the difference between the pI and your buffer pH increases.   |
|---------------------|--|
|                     | • In addition to pH, binding is also directly influenced by ionic strength (salt concentration). To ensure proper binding to resin, the starting sample salt concentration should be between 0-50mM for POROS 50 HQ and 0-150mM for POROS XQ.  |
| Flow rate           | • We recommend starting with a flow rate that allows for a 1-6 minute resident time (contact time) of a sample across the column. Flow rates can be increased or decreased significantly depending on sample concentration and complexity.   |
| Elution gradients   | • Running a salt gradient starting with your load salt concentration to 1M over 20 column volumes is recommended as a starting point. However, for some purifications >1M salt may be required to remove all bound components. Changes in pH can also be used. As an alternative to a salt gradient and pH gradient, a gradient combining shifts in pH and salt concentration could also be applied. |
|                     | • The slope at which a salt gradient is applied can have a significant impact on resolution. To optimize a purification protocol, use different gradient slopes throughout the protocol accordingly. Decrease the slope during points of interest to increase the resolution peaks and increase slopes at other times to minimize total purification time.   |
| Fraction collection | • The size of fractions to collect will be dependent on the complexity of your sample and the separation that can be achieved between sample components. This should be optimized for each purification. However, as a starting point, a fraction size of ~1/4–1/10 of the column volume can be used.  |
|                     | Pierce Biotechnology PO Box 117 (815) 968-0747 Thermofisher.com  |



#### **Standard purification protocol**

- 1. Equilibrate the column and all buffers to working temperature. Purifications can be performed at 2-30°C. Ensure that all solutions are degassed.
- 2. Prepare the purification system by washing pumps and filling tubing with buffer.
- 3. Connect the column to the system.
- 4. Equilibrate the column with 5-10 column volumes of the equilibration buffer.
- 5. Apply sample to the column. Under typical bind and elute protocols, the total protein load should not exceed 80% of the dynamic binding capacity of the column.
- 6. Wash the resin with >3 column volumes of wash buffer until the absorbance approaches baseline.
- 7. Elute bound molecules by applying a gradient from 0-100% elution buffer.
- 8. Maintain flow at 1M salt for 2-5 column volumes.
- 9. Wash the column with 3 column volumes of equilibration buffer.
- 10. Proceed to the "Cleaning-in-place and resin storage" section.

# **Purification: Flow-through**

A flow-through protocol is performed as a first-stage purification to quickly remove a large percentage of undesired proteins or as a final purification step in which your impurities are differentiated from your protein of interest, i.e. removal of HCP, DNA or virus from an antibody purification. In a flow-through protocol, the sample and equilibration buffer are adjusted to conditions where contaminant molecules will still bind to the resin, but your protein of interest will not. This is achieved by increasing the salt concentration and/or increasing the pH of the buffers to a point below the pI of your molecule of interest.

| Buffer composition | • When choosing buffer systems, consider molecule stability, binding optimization, and the ability of the buffer to control pH in the desired operating range (typically pH 7-9).  |
|--------------------|--|
| Conductivity       | • Typically, dynamic binding capacity decreases as load conductivity increases. POROS XQ and 50 HQ resins are salt-tolerant, allowing high-impurity dynamic binding capacity under higher conductivity load conditions. This reduces the need to dilute or buffer exchange column loads. The load conductivity should be 5-15mS/cm or about 50-150mM NaCl; however, the optimum buffer condition depends on the target molecule and buffer pH. |
| Capacity           | • The dynamic binding capacity must be experimentally determined for each impurity by using breakthrough analysis under the desired load pH and conductivity conditions.   |

- 1. Equilibrate the column and all buffers to working temperature. Perform purifications at 2-30°C. Ensure that all solutions are degassed.
- 2. Prepare the purification system by washing pumps and filling tubing with buffer.
- 3. Equilibrate the column with 5-10 column volumes of the equilibration buffer.
- 4. Begin applying sample while immediately collecting samples.

**Note:** The volume of sample that can be loaded will have to be determined experimentally as it is often difficult to determine when the dynamic binding capacity of the column will be exceeded by the unwanted binding materials.

- 5. Once sample has been fully applied, continue washing column with Equilibration Buffer until absorption reading returns to baseline.
- 6. Apply a gradient of 1M NaCl over 5 column volumes then hold at 1M NaCl for 2 column volumes or until material has been fully removed from column.
- 7. Wash the column with 3 column volume of equilibration buffer.
- 8. Proceed to the "Cleaning-in-place and resin storage" section.



# Cleaning-in-place and resin storage

Contaminants/impurities can often bind very tightly to POROS AEX resins in both bind/elute and flowthrough mode, so it is important to clean the POROS AEX resin sufficiently to increase column lifetime. The frequency at which this cleaning procedure is performed will be dependent on your sample, column performance, and regulatory needs. At a minimum perform the following clean-in-place procedure any time an increase in backpressure or decrease in performance is observed.

- 1. Wash the resin with 5 column volumes of 1M NaCl and 1M NaOH.
- 2. Wash the resin with 5 column volumes of ultrapure water.
- 3. Wash the resin with 5 column volumes of 1M acetic acid.
- 4. Wash the resin with 5 column volumes of ultrapure water.
- 5. Store resin in ultrapure water containing 20% ethanol at 2-30°C.

Note: If using the resin in flowthrough applications, additional cleaning may be needed.

# Troubleshooting

| Problem             | Cause                                  | Solution  |
|---------------------|--|---|
| Low protein yield   | Poor expression of soluble protein     | Optimize expression conditions. Possible conditions to<br>optimize are temperature of induction, length of induction,<br>media, and cell type |
|                     | Insufficient cell lysis and extraction | Optimize cell lysis protocol  |
|                     | Protein does not bind to the column    | Adjust binding buffer, modify salt concentration, or modify<br>pH to be further from pI of the protein of interest                            |
|                     | Flow rate was too fast                 | Decrease flow rate during binding to allow for greater residence time and increased binding   |
| Poor protein purity | Co-elution of contaminating            | Modify elution conditions or gradient   |
|                     | molecules                              | Use alternative IEX resin such as POROS HS or XS  |
|                     | Column was overloaded                  | Apply less protein extract onto the column  |
|                     | Column was dirty                       | Follow clean-in-place procedure to remove non-specifically bound impurities   |
| Slow column flow    | Extract was too viscous or highly      | Dilute lysate with equilibration buffer to decrease viscosity,  |
|                     | particulate                            | centrifuge lysate at higher speed to remove particulate, or   |
|                     |  | use Pierce <sup>TM</sup> Universal Nuclease for Cell Lysis (Product   |
|                     |  | No. 88/00) to eliminate DNA/RNA   |
| TT' - 1, 1, 1,      | Reprid switch in buffer tune           | Avoid rapid guitabas between organic and equeous  |
| High backpressure   | Kapid switch in burlet type            | solutions. Switch under reduced flow rates or over a  |
|                     |  | gradient  |
|                     | Presence of any amount of ethanol in   | Fully remove ethanol before packing or injecting sample   |
|                     | the slurry or column                   |   |
|                     | Compromised flow path                  | Characterize the pressure of the entire chromatography  |
|                     |  | system with 1) no column in place, 2) an empty column   |
|                     |  | with the column outlet plumbed to the waste, and 3) an  |
|                     |  | empty column with the column outlet plumbed back to the   |
|                     |  | chromatography system   |
|                     |  | Ensure the entire flow path is clear  |
|                     |  | Change in-line filters  |
|                     | Clogged frits                          | Change or clean frits   |
|                     |  | Run the column in reverse for 3 column volumes  |
|                     | Particle size gradient in the column   | Do not gravity settle POROS resin in the column before  |
|                     | caused by gravity settling             | Parask solumn   |
|                     | Resin was frozen                       | Do not freeze   |
|                     | Kesin was nozen                        |   |

Continued on next page

Pierce Biotechnology 3747 N. Meridian Road



| Continued from previo   | pus page   |  |
|---|--|--|
| Turbid column<br>effluent after   | Column frits are too large for the resin   | Use frits with <25µm pore size   |
| >3 column<br>volumes during<br>packing  | Defective O-ring or improperly<br>assembled column   | Disassemble column and inspect all components  |
| Column  | Column was underpacked   | Pack at higher flow rate/pressure  |
| qualification –<br>high asymmetry   | Salt injection was not optimized   | Verify the desired amount of salt is loaded by checking<br>the peak height and width   |
|   | System plumbing allowed for dilution of salt plug  | Characterize the salt plug through the chromatography<br>system at the qualification flow rate to understand how<br>the plug moves through the system with no packed<br>column in-line |
|   | Column needed more post-pack<br>conditioning to stabilize the packed<br>bed  | Equilibrate the column with 3-5 column volumes of packing solution in downflow at the operating flow rate  |
|   | Improper slug solution used  | Use the recommended column qualification conditions  |
| Column<br>qualification – low<br>asymmetry  | Water was used as the mobile phase   | Add some salt to the mobile phase to reduce the charge interaction between the salt and the resin  |
|   | Column was overpacked or packed inconsistently   | Repack the column following the recommended procedure  |
|   | Column not equilibrated long enough<br>with NaCl before salt injection   | Equilibrate with >4 column volumes if the packing solution is different from the qualification mobile phase  |
| Decreased<br>performance<br>• Increased band<br>spreading<br>• Dynamic binding<br>capacity decrease                             | Column fouling occurred due to<br>product precipitation or impurity,<br>irreversible binding of lipid material,<br>or other impurities | Clean the column   |
|   | Gradient was not correct   | Use a shallower gradient   |
|   |  | Lower flow rate  |
| <ul> <li>Recovery issues</li> </ul>   | Wrong type of gradient used  | Perform a pH gradient instead of salt gradient   |
| <ul> <li>Increased<br/>pressure drop</li> <li>Ghost peaks<br/>during other steps</li> <li>Protein<br/>separation pot</li> </ul> |  | Perform a salt gradient instead of a pH gradient   |
| achieved  |  |  |

8



| Related Thermo Scientific Products                                     |  |  |
|--|--|--|
| <b>POROS 50 HQ</b> , 1L (see thermofisher.com for larger pack sizes)   |  |  |
| POROS XQ, 1L   |  |  |
| POROS XS   |  |  |
| POROS PI 50µm, 1L  |  |  |
| POROS D 50µm, 1L   |  |  |
| B-PER <sup>™</sup> Bacterial Protein Extraction Reagent with Enzymes   |  |  |
| B-PER <sup>TM</sup> Bacterial Protein Extraction Reagent, 250mL        |  |  |
| <b>B-PER Bacterial Protein Extraction Reagent</b> , 500mL              |  |  |
| Halt <sup>TM</sup> Protease Inhibitor Cocktail (100X)                  |  |  |
| Pierce <sup>™</sup> BCA Protein Assay Kit, 1L                          |  |  |
| Zeba <sup>TM</sup> Desalting Chromatography Cartridges, $5 \times 5mL$ |  |  |
| Zeba <sup>TM</sup> Spin Desalting Columns, 5 × 10mL                    |  |  |
| Slide-A-Lyzer <sup>TM</sup> G2 Dialysis Cassettes                      |  |  |
|  |  |  |

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NÓ OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at thermofisher.com. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2016 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.

9