

# POROS and CaptureSelect products

## Your branding questions answered

In order to provide greater clarity to our customers, we recently retired the Life Technologies brand name and logo for what are now Thermo Scientific™ POROS™ and CaptureSelect™ products. We've kept a keen focus on ensuring that this simplification of our branding architecture will not impact the high level of service and support you've come to expect.

### **As part of our brand integration process, we're making the following changes:**

Newly manufactured products will be sold and shipped with updated branding and commercial documentation. This brand change won't impact the form, fit, or function of POROS\* and CaptureSelect\*\* products; it also won't affect the way you conduct business with us.

Expect to see the following branding changes:

- POROS and CaptureSelect product packaging, labels, and supporting documentation will now have the Thermo Scientific™ brand; however, there's no change to the products themselves
- Branding changes won't impact the ordering process, the Cat. Nos., or the way you do business with us

### **What is not changing:**

- Our products
- Our Cat. Nos.
- Our manufacturing
- Our legal entity
- Your order process
- Your purchase orders
- Your invoice
- Your pricing
- Your packing slips
- Your shipping and delivery times

There's no change to how you do business with us as your vendor—we'll continue serving your bioproduction needs.

### **Where you'll see new branding:**

- Product packaging
- Shipping boxes
- Documentation and literature
- Certificates of analysis
- Material safety data sheets
- Marketing material and advertising

# You still get the same products and the same great features

## **What's happening to the Life Technologies brand?**

We have retired the Life Technologies brand. CaptureSelect products now have the Thermo Scientific brand.

## **What's happening to the Applied Biosystems™ brand?**

We didn't retire the brand; however, POROS purification products now have the Thermo Scientific brand.

## **What does this change mean to me?**

The same high-quality products are still developed, manufactured, and supported by the same life science team. There are no changes to our products.

## **When will I see these changes? When will the changes be complete?**

We've already begun shipping products with the Thermo Scientific brand. Over the next year you may see a combination of existing and new packaging as we exhaust our preexisting inventory to minimize impact on the environment.

## **Why are these changes being made?**

We're simplifying our brand architecture to create greater clarity for our customers. We believe our value to you lies in the well-respected brands you know and trust, as well as the choice afforded to you by one of the most innovative life science portfolios in the industry. That's not changing.

## **Where can I learn more about these changes?**

Go to [thermofisher.com/purificationbrandchange](https://thermofisher.com/purificationbrandchange) for more detailed information on the brand change, or [thermofisher.com/bioproduction](https://thermofisher.com/bioproduction) to learn more about all of our bioproduction products.

## **How should I reference your company in my publications?**

Please refer to the company as Thermo Fisher Scientific. When listing products in materials and methods sections, please use the Thermo Scientific brand before the product name.

## **Are there brand changes to other products?**

The brand and logo changes only relate to POROS and CaptureSelect products.

## **Does my vendor change?**

No. There are no changes to your vendor as a result of this rebranding.

## **Will transactional documents change?**

Invoices and other transactional documents will have the logo of Thermo Fisher Scientific, but will reference the appropriate legal entity. This brand transition doesn't affect our legal entities.

## Product information

### **Are catalog numbers changing?**

No. Our Cat. Nos. are not changing as a result of this rebranding.

### **Are there changes to product formulations?**

No. Our manufacturing processes are not changing as a result of the brand simplification.

### **Are there changes in quality control or testing?**

No. Our quality processes are not changing as a result of the brand simplification.

### **Are there changes to product labels?**

Yes. The brand name and logo on some of our product labels will change. However, Life Technologies will remain as the legal entity/manufacturer of the product. There are no changes to our products.

### **Are there changes to product packaging?**

Yes. The brand name and logo on some of our product packaging will change. We're depleting our supply of existing brand packaging to minimize waste, so over the next year you may see a combination of existing and new packaging.

### **Are there changes to certificates of analysis?**

Yes. The brand name and logo on some of our certificates will change. However, these documents will retain Life Technologies as the legal entity/manufacturer, as appropriate.

### **Are there changes to product literature?**

Yes. Product literature will have an updated Thermo Scientific brand.

### **Are there changes to shipping boxes?**

Yes. The brand name and logo on our shipping boxes will have the Thermo Fisher Scientific company name. We're depleting our supply of existing shipping boxes to minimize waste, so you may see a combination of existing and new shipping boxes.

## Doing business with us

### **Should I change the name of the vendor in my purchasing systems?**

No. There are no changes to our legal entity names or contact information; please continue to submit purchase orders and payments to the same vendor you do today.

### **Are there changes to ordering processes?**

No. There are no changes to ordering as a result of this rebranding.

### **Should I change the name of the vendor/legal entity on my order?**

No. Please continue to use the existing vendor/legal entity name you do today.

### **Are there changes to my existing contract?**

No. There are no changes to your existing contract as a result of the brand change.

### **Are there changes to the terms and conditions?**

No. There are no changes to the terms and conditions as a result of this rebranding.

### **Are there changes to my pricing?**

No. There are no changes to pricing as a result of this rebranding.

### **Are there changes to freight or handling?**

No. There are no changes to freight or handling as a result of this rebranding.

### **Are there changes to my authorized distributor?**

No. Please continue to order as usual.

Find out more at  
[thermofisher.com/purificationbrandchange](http://thermofisher.com/purificationbrandchange)

**ThermoFisher**  
SCIENTIFIC

\* POROS products are a **Pharmaceutical Grade Reagent. For Manufacturing and Laboratory Use Only.** Some POROS products <1 L may be for **Research Use Only. Not for use in diagnostic procedures.** See individual product labels for details.

\*\* CaptureSelect products are **For In-Process Quality Control testing.**