

Changes to Gibco product use statement, logo, and packaging

FAQs

What's happening to the Gibco™ brand logo?

The old logo includes a tagline that reads “Gibco by Life Technologies.” In November 2017, we will be removing this tagline, leaving just the Gibco brand logo.

How does this change affect me?

There will be no changes to our products, nor how they're developed, manufactured, and supported. The only changes will take place on our documentation, packaging, and labeling, including the product use statement on some products. Depending upon your processes, this may require you to change your specifications.

When will I see these changes?

You should begin to see changes in November 2017. You may see a combination of existing and new packaging as we exhaust our preexisting inventory.

Why are these changes being made?

Based on customer feedback, we're clarifying the logo and use statements. We're not changing the products nor how you conduct business with us.

Where can I learn more about these changes?

Go to thermofisher.com/gibcolabelchanges for more detailed information on the product label statement and branding changes, or thermofisher.com/bioproduction to find out more about our bioproduction products.

How should I reference your company in my publications?

Please refer to the company as Thermo Fisher Scientific.

Are there label statement or brand changes to other products?

The label statement and logo changes only relate to Gibco™ products.

Does my vendor change?

No. There will be no changes to our vendor name nor the legal entity.

Will transactional documents change?

Invoices and other transactional documents will continue to reference the appropriate legal entity. The label statement and brand change doesn't affect our legal entities.

Who should I contact if I have questions?

Please email techsupport@thermofisher.com or call **800-955-6288** opt 5, ext. 40562, in US/Canada or **00-800-5345-5345** option 3 in Europe. For all other regions, please refer to our website for your local phone number.

Find out more at thermofisher.com/gibcolabelchanges

ThermoFisher
SCIENTIFIC