



September 2025

Change Notification Letter

Dear valued customer,

QIAGEN is dedicated to meeting the needs of our customers, and we continually strive to provide the best value in products and services. As such, we are writing to inform you of a change impacting the following products.

Affected Products

<u>Catalog No.</u>	<u>Product</u>
31014	Ni-NTA Spin Column (50)
31314	Ni-NTA Spin Kit (50)

Description of Change

We exchange the color of the Ni-NTA Spin column from green color to transparent. The color has no impact on product performance in terms of quality or quantity of the isolated protein. The kit might include green or transparent Ni-NTA Spin columns within the transition phase.

This change has no impact on the fit, form, or functionality of the product.

If you would like more information about this change notification, please contact QIAGEN Technical Services at support.qiagen.com

Sincerely,

Your QIAGEN Quality Team

Trademarks: QIAGEN®, Sample to Insight® (QIAGEN Group). Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are not to be considered unprotected by law.

1138636 HB-3753-001 © 2025 QIAGEN, all rights reserved.