

Important customer notification

Pub. No. MAN0019724 Rev. A.0

The purification resin in the Antibody and the Protein Labeling kits has been changed to our new Zeba™ Dye and Biotin Removal Column. The Zeba™ Dye and Biotin Removal Column offers superior purification efficiency and ease compared to our previous resin.

You can find the user guide for the new purification system at **thermofisher.com**.

You can still purchase the previous resin and purification system for the Antibody and the Protein Labeling kits as a separate kit as listed below:

Antibody Conjugate Purification Kit for 20–50 µg (Cat. No. A33087)

Antibody Conjugate Purification Kit for 50–100 µg (Cat. No. A33088)

Antibody Conjugate Purification Kit for 0.5–1 mg (Cat. No. A33086)

For Research Use Only. Not for use in diagnostic procedures.

ThermoFisher
SCIENTIFIC



Life Technologies Corporation | 29851 Willow Creek | Eugene, OR 97402

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit thermofisher.com/support
thermofisher.com

ThermoFisher
SCIENTIFIC