



ChromaLINK[®] Biotin Protein Labeling Kit

B-9007-105

[Product Images](#)



Short Description

The ChromaLINK Biotin Protein Labeling Kit includes everything you need to biotinylate your protein: buffers, reagents, spin columns, and an easy-to-follow protocol and online calculator. With the ChromaLINK Biotin Protein Labeling Kit you can biotinylate and purify up to 5 antibodies or other proteins ranging in molecular weight from 20–200 kDa and from as little as 25 µg to as much as 1 mg.

Zeba and Thermo Scientific are trademarks of Thermo Fisher Scientific Inc.

Additional Information

Conjugation Targets	Antibody
Label/Modifier Type	biotin
Reactivity	Amine
Recommended Storage	2° - 8°C — Do Not Freeze
Applications	Antibody Labeling

Products in this set

ChromaLINK® Biotin Protein Labeling Kit

TriLink's large scale ChromaLink® One-Shot Antibody Biotinylation Kit provides convenient, consistent and measurable biotinylation of up to 2 mg of antibody. The kit contains ChromaLink® Biotin, which incorporates a UV-traceable chromophore based on Chromalink® technology to enable reproducibility in your biotinylation process. Now you can measure the degree of biotinylation in minutes, not hours, without the standard curves required for HABA/avidin and fluoro-reporter assays. The One-Shot kit contains everything needed to biotinylate your antibody: buffers, reagents, spin columns, and an easy-to-follow protocol and online MSR calculator.



