



# ChromaLINK® One-Shot™ Antibody Biotinylation Kit

## B-9007-009

[Product Images](#)



## Short Description

---

The ChromaLINK One-Shot Antibody Biotinylation Kit provides convenient, consistent, and measurable biotinylation of 100 µg of antibody. Each kit contains ChromaLINK Biotin, which incorporates a novel UV-traceable chromophore in the linker arm 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 the HABA/Avidin and Fluoro-Reporter assays. The ChromaLINK One-Shot kit contains everything needed to biotinylate your antibody: buffers, reagents, desalting columns, and an easy-to-follow protocol and online biotin incorporation calculator.

## Additional Information

---

Conjugation Targets	Antibody
Label/Modifier Type	biotin
Reactivity	Amine
Recommended Storage	2° - 8°C — Do Not Freeze
Applications	Antibody Labeling, In Situ Proximity Ligation
Unit Size	1 Kit

## Products in this set

---



**ChromaLINK® One-Shot™ Antibody Biotinylation Kit**

