



ChromaLINK® Digoxigenin One-Shot™ Antibody Labeling Kit

B-9014-009

[Product Images](#)



Short Description

The ChromaLINK Digoxigenin One-Shot Antibody Labeling Kit provides convenient, consistent, and measurable digoxigenin labeling of 100 µg of antibody. Each kit contains ChromaLINK Digoxigenin, which incorporates a novel UV-traceable chromophore in the linker arm to enable reproducibility in your antibody labeling process. With a simple, non-destructive UV scan you can now quantify digoxigenin incorporation to ensure reproducible incorporation of the optimal number of haptens per antibody. Each ChromaLINK One-Shot kit contains everything needed to label your antibody: buffers, reagents, desalting columns, and an easy-to-follow protocol and online DIG incorporation calculator.

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

Additional Information

Conjugation Targets	Antibody
Label/Modifier Type	Digoxigenin
Reactivity	Amine
Recommended Storage	2° - 8°C – Do Not Freeze
Applications	In situ hybridization, Antibody Labeling, In Situ Proximity Ligation

Products in this set



ChromaLINK® Digoxigenin One-Shot™ Antibody Labeling Kit

