

154Sm Metal Conjugation Kit - Lightning-Link® ab269837

1 Image

Overview

Product name 154Sm Metal Conjugation Kit - Lightning-Link®

Product overview 154Sm Metal Conjugation Kit - Lightning-Link® (ab269837) provides an easy-to-use, one step procedure that allows researchers to label proteins, peptides and other biomolecules containing primary amines with 154Sm with only 30 seconds hands-on time; furthermore conjugates are ready to use in less than twenty minutes.

This product is manufactured by Expedeon, an Abcam company. Expedeon product code M154-0010 was previously called Lightning-Link 154Sm, 1 x 10 µg and is the same as the 10 µg size of this product. Expedeon product code M154-0100 was previously called Lightning-Link 154Sm, 1 x 100 µg and is the same as the 100 µg size.

The antibody to be labeled should be purified, in an appropriate buffer for conjugation and at a suitable concentration.

Learn about buffer compatibility below; for incompatible buffers and low antibody concentrations, use our rapid [antibody purification and concentration kits](#). Use the [FAQ](#) to learn more about the technology, or about conjugating other proteins and peptides to 154Sm.

Notes

Buffer Requirements for Conjugation

Buffer should be pH 6.5-8.5.

Compatible buffer constituents

If a concentration is shown, then the constituent should be no more than the concentration shown. If several constituents are close to the limit of acceptable concentration, then this can inhibit conjugation.

50mM / 0.6% Tris ¹	0.1% BSA ²	50% glycerol
0.1% sodium azide	PBS	Potassium phosphate
Sodium chloride	HEPES	Sucrose
Sodium citrate	EDTA	Trehalose

¹ Tris buffered saline is almost always ≤ 50 mM / 0.6%

² BSA can also interfere with the use of the conjugated antibody in tissue staining.

Incompatible buffer constituents

Thiomerosal	Proclin	Glycine
Arginine	Glutathione	DTT

If a constituent of the buffer containing your antibody or protein is not listed above, please check the [FAQ](#) or [contact us](#).

Only purified antibodies are suitable for use, ie. where other proteins, peptides, or amino acids are not present: antibodies in ascites fluid, serum or hybridoma culture media are incompatible.

Storing and handling conjugation kits

Lyophilized Lightning-Link[®] components are hygroscopic.

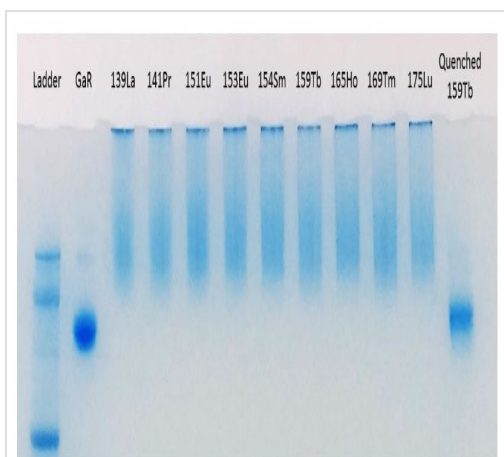
Kits are intentionally shipped at ambient temperature with silica gel to avoid exposure to moisture. Upon receipt, store the kit frozen and protect from moisture. Before opening the outer container, allow the lyophilized components to reach room temperature to minimize condensation.

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	100 µg
ab274207 - 154Sm mix (Lyophilized)	1 x 100µg
ab273994 - Modifier reagent	1 x 200µl
ab274133 - Quencher reagent	1 x 200µl

Images



Gel Image Showing Lightning-Link® Metal Labeling Kit Antibody Labeling Efficiency.

Goat Anti-Rabbit was labeled with a selection of Lightning-Link® metal labels at 10 µg single reaction vials. 2 µg conjugates per well were run on a 4-12% Bis-Tris gel, under non-reducing conditions for 20 minutes at 50V and 1 hour at 120V in 1x MOPS. The gel was stained with InstantBlue™ for 1 hour on the gel rocker. The gel showed no free antibody post conjugation for all conjugates except when the reaction was quenched at t=0, where no conjugation was observed.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors