

Atto 565 Conjugation Kit (Fast) - Lightning-Link® ab269897

[3 References](#) [1 Image](#)

Overview

Product name Atto 565 Conjugation Kit (Fast) - Lightning-Link®

Product overview Atto 565 Conjugation Kit (Fast) - Lightning-Link® (ab269897) provides an easy-to-use, one step procedure that allows researchers to covalently label proteins, peptides and other biomolecules containing primary amines with Atto 565 with only 30 seconds hands-on time; furthermore conjugates are ready to use in less than twenty minutes.

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 Atto 565.

Custom size conjugation kits up to 100 mg are available on demand. Please contact us to discuss your requirements.

Notes This product is manufactured by Expedeon, an Abcam company, and was previously called Lightning-Link® Rapid Atto 565 Labeling Kit. 351-0005 is the same as the 100 µg size. 351-0010 is the same as the 3 x 100 µg size. 351-0030 is the same as the 3 x 10 µg size.

Amount and volume of antibody for conjugation to Atto 565

<i>Kit size</i>	<i>Recommended amount of antibody¹</i>	<i>Maximum amount of antibody</i>	<i>Maximum antibody volume²</i>
3 x 10 µg	3 x 10 µg	3 x 20 µg	3 x 10 µL
100 µg	100 µg	200 µg	100 µL
3 x 100 µg	3 x 100 µg	3 x 200 µg	3 x 100 µL

¹ Using the maximum amount of antibody may result in less labeling per antibody.

² Ideal antibody concentration is 1mg/ml. 0.5 - 1 mg/ml can be used if the maximum antibody volume is not exceeded. Antibodies > 2 mg/ml or < 0.5 mg/ml should be diluted /concentrated.

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.

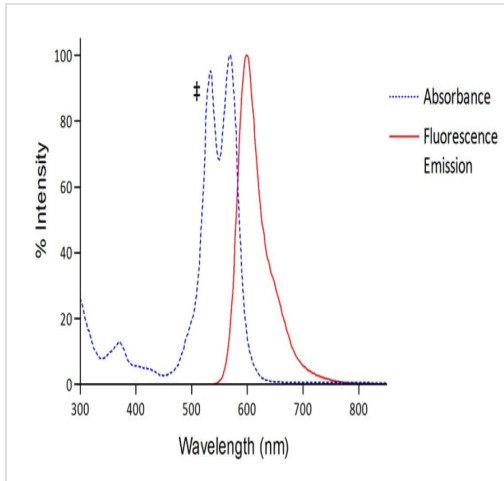
Properties

Storage instructions

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

Components	100 µg	3 x 10 µg	3 x 100 µg
ab274068 - Atto 565 mix (Lyophilized)	1 x 100µg	3 x 10µg	3 x 100µg
ab273994 - Modifier reagent	1 x 200µl	1 x 200µl	1 x 200µl
ab273995 - Quencher reagent	1 x 200µl	1 x 200µl	1 x 200µl

Images



†† This peak is an absorbance peak rather than a fluorescence excitation peak, so the dye will not fluoresce if excitation wavelengths in this range are used.

Atto 565 Conjugation Kit (Fast) - Lightning-Link®
absorbance emission graph

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

Our Promise 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