

Atto 633 Conjugation Kit (Fast) - Lightning-Link® ab269898

[7 References](#) [2 图像](#)

概述

产品名称

Atto 633偶联试剂盒(Fast) - Lightning-Link®

产品概述

Atto 633 Conjugation Kit (Fast) - Lightning-Link® (ab269898) provides an easy-to-use, one step procedure that allows researchers to covalently label proteins, peptides and other biomolecules containing primary amines with Atto 633 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.

The excitation and emission wavelengths for Atto 633 are Ex: 630nm, Em: 651nm.

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 633.

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

说明

This product is manufactured by Expedeon, an Abcam company, and was previously called Lightning-Link® Rapid Atto 633 Labeling Kit. 353-0015 is the same as the 1 mg size. 353-0010 is the same as the 3 x 100 µg size. 353-0030 is the same as the 3 x 10 µg size. 353-0005 is the same as the 100 µg size.

Amount and volume of antibody for conjugation to Atto 633

<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	1 x 100 µg	1 x 200 µg	1 x 100 µL
3 x 100 µg	3 x 100 µg	3 x 200 µg	3 x 100 µL
1 mg	1 x 1 mg	1 x 2 mg	1 x 1 mL

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 > 2mg/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.

性能

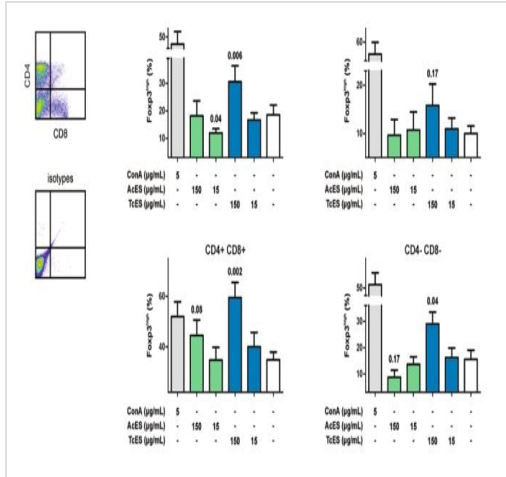
存放说明

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

组件	100 µg	1 mg	3 x 10 µg	3 x 100 µg
ab274070 - Atto 633 mix (Lyophilized)	1 x 100µg	1 x 1mg	3 x 10µg	3 x 100µg
ab273994 - Modifier reagent	1 x 200µl	1 x 200µl	1 x 200µl	1 x 200µl

组件	100 µg	1 mg	3 x 10 µg	3 x 100 µg
ab273995 - Quencher reagent	1 x 200µl	1 x 200µl	1 x 200µl	1 x 200µl

图片

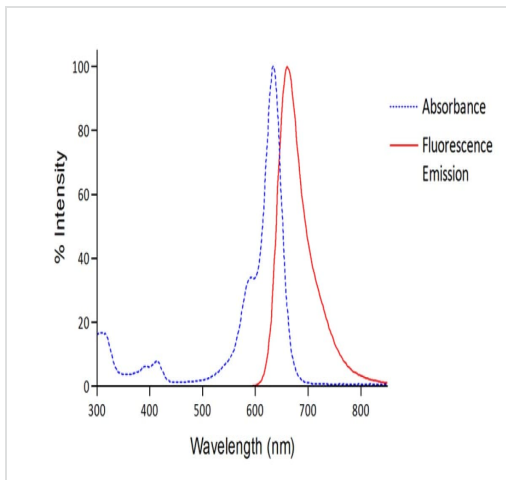


Flow Cytometry - Atto 633 Conjugation Kit (Fast) - Lightning-Link (ab269898)

Image from Junginger, Johannes et al., Scientific reports vol. 7, 1 10310. 4 Sep. 2017, doi:10.1038/s41598-017-10677-4. Reproduced under the Creative Commons license <https://creativecommons.org/licenses/by/4.0/>

Junginger, Johannes et al used Atto 633 Conjugation Kit - Lightning-Link® (ab269898) as part of examining zoonotic intestinal helminths. They used the kit to conjugate Atto 633 to anti-canine CD4 antibody for use in flow cytometry.

Three-colour flow cytometry revealed the TcES-associated increase in Fopx3high lymphocytes to be associated with CD4+, CD4+ CD8+ double-positive and CD4- CD8- double-negative subsets, while this effect was lower in CD8+ T cells. Compared to TcES, treatment with AcES at 150 µg/mL was associated with a much lower elevation in Fopx3high expression by lymphocytes and CD4+ CD8+ T cells.



Atto 633 Conjugation Kit (Fast) - Lightning-Link® absorbance emission graph

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.cn/abpromise> or contact our technical team.

Terms and conditions

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