

# Alexa Fluor® 594 Conjugation Kit (Fast) - Lightning-Link® ab269822

## 5 References

### Overview

**Product name** Alexa Fluor® 594 Conjugation Kit (Fast) - Lightning-Link®

**Product overview** Alexa Fluor® 594 Conjugation Kit / Alexa Fluor® 594 Labeling Kit (ab269822) uses a simple and quick process for Alexa Fluor® 594 labeling / conjugation of antibodies. It can also be used to conjugate other proteins or peptides. Learn about our [antibody labeling kits and their advantages](#).

To conjugate an antibody to Alexa Fluor® 594 using this kit:

- add modifier to antibody and incubate for 15 mins
- add quencher and incubate for 5 mins

The Alexa Fluor® 594 conjugated antibody can be used immediately in WB, ELISA, IHC etc. No further purification is required and 100% of the antibody is recovered for use.

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 Alexa Fluor® 594.

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 Alexa Fluor® 594 Labeling Kit. 335-0015 is the same as the 1 mg size. 335-0010 is the same as the 3 x 100 ug size. 335-0030 is the same as the 3 x 10 ug size. 335-0005 is the same as the 100 µg size.

### Amount and volume of antibody for conjugation to Alexa Fluor® 594

| <i>Kit size</i> | <i>Recommended amount of antibody<sup>1</sup></i> | <i>Maximum amount of antibody</i> | <i>Maximum antibody volume<sup>2</sup></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 mg | 1 x 1 mg | 1 x 2 mg | 1 x 1 mL |
|------|----------|----------|----------|

<sup>1</sup> Using the maximum amount of antibody may result in less labelling per antibody.

<sup>2</sup> 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 <sup>1</sup> | 0.1% BSA <sup>2</sup> | 50% glycerol        |
|-------------------------------|-----------------------|---------------------|
| 0.1% sodium azide             | PBS                   | Potassium phosphate |
| Sodium chloride               | HEPES                 | Sucrose             |
| Sodium citrate                | EDTA                  | Trehalose           |

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

<sup>2</sup> 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<sup>®</sup> 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.

Alexa Fluor<sup>®</sup> is a registered trademark of Molecular Probes, Inc, a Thermo Fisher Scientific Company. The Alexa Fluor<sup>®</sup> dye included in this product is provided under an intellectual property license from Life Technologies Corporation. As this product contains the Alexa Fluor<sup>®</sup> dye, the purchase of this product conveys to the buyer the non-transferable right to use the purchased product and components of the product only in research conducted by the buyer (whether the buyer is an academic or for-profit entity). As this product contains the Alexa Fluor<sup>®</sup> dye the sale of this product is expressly conditioned on the buyer not using the product or its components, or any materials made using the product or its components, in any activity to generate revenue, which may include, but is not limited to use of the product or its components: in manufacturing; (ii) to

provide a service, information, or data in return for payment (iii) for therapeutic, diagnostic or prophylactic purposes; or (iv) for resale, regardless of whether they are sold for use in research. For information on purchasing a license to this product for purposes other than research, contact Life Technologies Corporation, 5781 Van Allen Way, Carlsbad, CA 92008 USA or [outlicensing@thermofisher.com](mailto:outlicensing@thermofisher.com).

## Properties

---

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

| Components                  | 100 µg    | 1 mg      | 3 x 10 µg | 3 x 100 µg |
|-----------------------------|-----------|-----------|-----------|------------|
| ab274046 - Alexa Fluor 594  | 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  |
| ab273995 - Quencher reagent | 1 x 200µl | 1 x 200µl | 1 x 200µl | 1 x 200µl  |

**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