

Texas Red® Conjugation Kit (Fast) - Lightning-Link® ab195225

7 References [画像数 3](#)

製品の概要

製品名 Texas Red® Conjugation Kit (Fast) - Lightning-Link®

製品の概要 Texas Red® Conjugation Kit / Texas Red Labeling Kit ab195225 uses a simple and quick process for Texas Red® 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 Texas Red® using this kit:

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

The Texas Red 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 Texas Red®.

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

特記事項

This product is manufactured by Expedeon, an Abcam company, and was previously called Lightning-Link® Rapid Texas Red® Labeling Kit. 315-0005 is the same as the 100 ug size. 315-0010 is the same as the 3 x 100 ug size. 315-0030 is the same as the 3 x 10 ug size. 315-0015 is the same as the 1 mg size.

Amount and volume of antibody for conjugation to Texas Red®

<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 labelling 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.

製品の特性

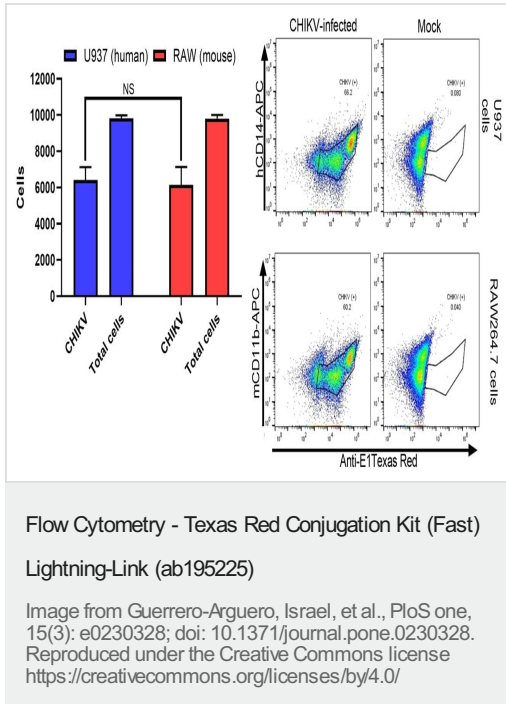
保存方法

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

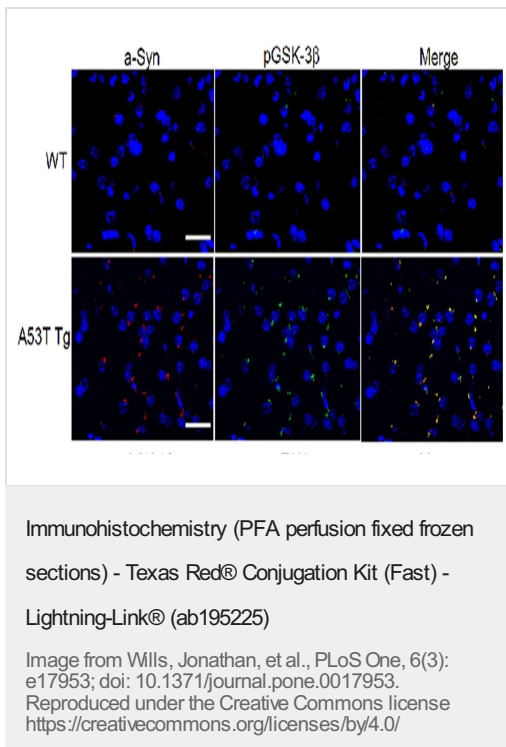
内容	1 mg	100 µg	3 x 10 µg	3 x 100 µg
ab273994 - Modifier reagent	1 x 200µl	1 x 200µl	1 x 200µl	1 x 200µl
ab273995 - Quencher reagent	1 x 200µl	1 x 200µl	1 x 200µl	1 x 200µl

内容	1 mg	100 µg	3 x 10 µg	3 x 100 µg
ab274004 - Texas Red Conjugation Mix	1 x 1mg	1 x 100µg	3 x 10µg	3 x 100µg

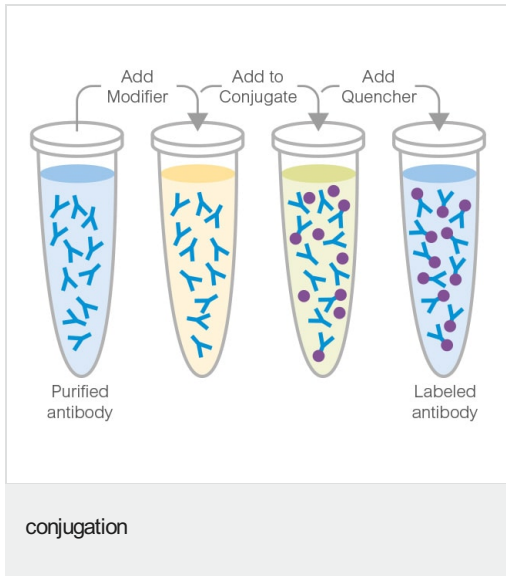
画像



Guerrero-Arguero, Israel, et al used Texas Red® Conjugation Kit (Fast) - Lightning-Link® (ab195225) as part of examining the rate of productive Chikungunya virus (CHIKV) replication via flow cytometry. They used the kit to conjugate Texas Red® to anti-Chikungunya E1 glycoprotein antibody for use in flow cytometry. PMA-differentiated U937 and RAW264.7 cells display CHIKV envelope proteins at 6 hpi. PMA = phorbol 12-myristate 13-acetate PMA-differentiated U937 and RAW264.7 macrophages were exposed for 2 hours to CHIKV and then fixed and assayed at 8 hpi using flow cytometry and an anti-E1 protein fluorophore-conjugated monoclonal antibody. Data show mean values of three independent experiments with a total of n = 9, MOI = 1. Statistical significance was determined using multiple t-test corrected using Holm-Sidak method. *P<0.05; NS, not significant.



Wills, Jonathan, et al used Texas Red® Conjugation Kit (Fast) - Lightning-Link® (ab195225) as part of examining the aggregation and co-localization of α-Syn with pGSK-3? and p-Tau in the striatum of A53T α-Syn mutant mice (A53T Tg) and wild-type mice (WT). They used the kit to conjugate Texas Red® to anti-α-Syn antibody for use in immunohistochemistry (PFA perfusion fixed frozen sections). Right panels constitute merged image of left panels. Sections of striatum of A53T Tg and age-matched WT mice were stained with anti-α-Syn antibody conjugated to Texas Red (red) and anti-p-GSK3β conjugated to FITC (green). Nuclei were stained with DAPI (blue).



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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