

# ReadiLink™ Rapid XFD555 Antibody Labeling Kit \*XFD555 Same Structure to Alexa Fluor™ 555\*

Catalog number: 1276 Unit size: 2x50 ug Labelings

Component	Storage	Amount
Component A: XFD555	Freeze (< -15 °C), Minimize light exposure	2 vials (One vial is for 50 μg protein)
Component B: Reaction Buffer	Freeze (< -15 °C), Minimize light exposure	1 vial (20 μL)
Component C: TQ™-Dyed Quench Buffer	Freeze (< -15 °C), Minimize light exposure	1 vial (20 μL)

#### **OVERVIEW**

XFD555 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 555 (Alexa Fluor® is the trademark of ThermoFisher). Readilink™ protein labeling technology is the most robust and convenient tool for preparing fluorescent antibody conjugates used for fluorescence imaging and flow cytometry applications. ReadiLink™ Rapid XFD555 Antibody Labeling Kit provides the most convenient tool for making XFD555-labeled antibody conjugates. XFD555 dye used in this ReadiLink™ kit is reasonably stable and shows good reactivity and selectivity with protein amino groups. The kit has all the essential components for labeling ~2x50 ug antibody. Each of the two vials of XFD555 dye provided in the kit is optimized for labeling ~50 µg antibody. ReadiLink™ Rapid XFD555 Antibody Labeling Kit requires minimal hands-on time. The prepared XFD555 conjugates (with the kit) are ready to use for fluorescence imaging and flow cytometry applications without further purifications needed. It provides the most convenient method to prepare XFD555-labeled antibodies.

## AT A GLANCE

#### Important

Warm all the components and centrifuge the vials briefly before opening, and immediately prepare the required solutions before starting your conjugation. The following protocol is for recommendation.

# PREPARATION OF WORKING SOLUTION

### Protein working solution (Solution A)

For labeling 50  $\mu g$  of protein (assuming the target protein concentration is 1 mg/mL), mix 5  $\mu L$  (10% of the total reaction volume) of Reaction Buffer (Component B) with 50  $\mu L$  of the target protein solution.

**Note** If you have a different protein concentration, adjust the protein volume accordingly to make  $\sim$ 50  $\mu$ g of protein available for your labeling reaction.

**Note** For labeling 100  $\mu$ g of protein (assuming the target protein concentration is 1 mg/mL), mix 10  $\mu$ L (10% of the total reaction volume) of Reaction Buffer (Component B) with 100  $\mu$ L of the target protein solution.

Note The protein should be dissolved in 1X phosphate buffered saline (PBS), pH 7.2 - 7.4; if the protein is dissolved in glycine buffer, it must be dialyzed against 1X PBS, pH 7.2 - 7.4, or use ReadiUse™ Bio-Gel P-6 spin column (cat# 60500 from AAT Bioquest) to remove free amines or ammonium salts (such as ammonium sulfate and ammonium acetate) that are widely used for protein precipitation.

**Note** Impure antibodies or antibodies stabilized with bovine serum albumin (BSA) or gelatin will not be labeled well.

Note For optimal labeling efficiency, a final protein concentration range of 1 - 2 mg/mL is recommended, with a significantly reduced conjugation efficiency at less than 1 mg/mL.

# SAMPLE EXPERIMENTAL PROTOCOL

### Run conjugation reaction

 Add the protein working solution (Solution A) to ONE vial of labeling dye (Component A-XFD555), and mix them well by repeatedly pipetting for a few times or vortex the vial for a few seconds.

**Note** If labeling 100  $\mu$ g of protein, use both vials (Component A) of labeling dye by dividing the 100  $\mu$ g of protein into 2 x 50  $\mu$ g of protein and reacting each 50  $\mu$ g of protein with one vial of labeling dye. Then combine both vials for the next step.

 Keep the conjugation reaction mixture at room temperature for 30 -60 minutes.

**Note** The conjugation reaction mixture can be rotated or shaken for longer time if desired.

## Stop Conjugation reaction

- Add 5 μL (for 50 μg protein) or 10 μL (for 100 μg protein) which is 10% of the total reaction volume of TQ™-Dyed Quench Buffer (Component C) into the conjugation reaction mixture; mix well.
- Incubate at room temperature for 10 minutes. The labeled protein (antibody) is now ready to use.

# Storage of Protein Conjugate

The protein conjugate should be stored at > 0.5 mg/mL in the presence of a carrier protein (e.g., 0.1% bovine serum albumin). For longer storage, the protein conjugates could be lyophilized or divided into single-used aliquots and stored at  $\leq -20^{\circ}$ C.

# DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.