

# ab312831 – APC-iFluor® 700 Conjugation Kit

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: [www.abcam.com/ab312831](http://www.abcam.com/ab312831)

## Storage and Stability:

Upon receipt, store the kit at 4 °C. When stored properly, the kit should be stable for six months. Alternatively, Buccutite™ MTA can be stored at -20 °C. Do not freeze Buccutite™ FOL-Activated APC-iFluor™ 700 , Reaction Buffer . Warm all the components and centrifuge the vials briefly before opening, and immediately prepare the required solutions before starting your conjugation. The following SOP is an example for labeling goat anti-mouse IgG antibody.

## Materials Supplied

Item	Quantity	Storage Condition
Buccutite™ FOL-Activated APC-iFluor™ 700	2 vials	-20°C
Buccutite™ MTA	1 vials	-20°C
Reaction Buffer	1 vials	-20°C

## Reagent Preparation

**Antibody working solution:** For labeling 25 µg antibody (assuming the target antibody concentration is 1 mg/mL), mix 1.25 µL (5% of the total reaction volume) of Reaction Buffer with 25 µL of the target antibody solution.

**Note:** If you have a different concentration, adjust the antibody volume accordingly to make ~25 µg antibody available for your labeling reaction.

**Note:** The antibody should be dissolved in 1X phosphate buffered saline (PBS), pH 7.2-7.4; If the antibody is dissolved in glycine buffer, it must be dialyzed against 1X PBS, pH 7.2-7.4, or use Amicon Ultra-0.5, Ultracel-10 Membrane, 10 kDa ( Cat. # UFC501008 from Millipore) to remove free amines or ammonium salts (such as ammonium sulfate and ammonium acetate) that are widely used for antibody precipitation.

**Note:** Impure antibodies or antibodies stabilized with bovine serum albumin (BSA) or gelatin will not be labeled well. For BSA removal, we recommend BSA Removal Kit (ab173231). For purification and concentration we recommend Antibody Concentration And Clean-Up Kit (ab102778).

**Note:** The antibody –Buccutite™ MTA reaction efficiency is significantly reduced if the antibody concentration is less than 1 mg/mL. For optimal labeling efficiency the final antibody concentration range of 1-10 mg/mL is recommended.

**Buccutite™ MTA working solution:** Add 10 µL DMSO (Not provided) into the vial of Buccutite™ MTA

**Buccutite™ FOL-Activated APC-iFluor™ 700 working solution:** Add 50 µL ddH<sub>2</sub>O into the vial of Buccutite™ FOL-Activated APC-iFluor™ 700

## Assay Procedure

### Run Antibody-Buccutite™ MTA reaction

1. Add 2.5 µL of Buccutite™ MTA working solution into antibody working solution, and mix them well by repeatedly pipetting for a few times or vortex the vial for a few seconds.
2. Keep the antibody- Buccutite™ MTA reaction mixture at room temperature for 30 - 60 minutes.

**Note:** The antibody-Buccutite™ MTA reaction mixture can be rotated or shaken for longer time if desired.

### Make antibody-APC-iFluor™ 700 conjugation

1. Add 50 µL of Buccutite™ FOL-Activated APC-iFluor™ 700 working solution with Antibody-Buccutite™ MTA solution, mix well by repeatedly pipetting for a few times or vortex the vial for a few seconds.
2. Incubate for 1 to 2 hours
3. The antibody-APC-iFluor™ 700 conjugate is now ready to use.

**Note:** For immediate use, the antibody-APC-iFluor™ 700 conjugate need be diluted with the buffer of your choice.

**Note:** For longer term storage, antibody-APC-iFluor™ 700 conjugate solution need be concentrated or freeze dried

### Storage of Antibody-APC-iFluor™ 700 Conjugate

The antibody conjugate should be stored in the presence of a carrier protein (e.g., 0.1% bovine serum albumin) and 0.02-0.05% sodium azide. The Ab-APC-iFluor™ 700 conjugate solution could be stored at 4 °C for two months without significant change and kept from light