

Goat Anti-Mouse IgG, F(ab')₂ Fragment, Fc Gamma-Specific, Highly Cross-Adsorbed, CF[®]640R Conjugate



Highly cross-adsorbed F(ab')₂ fragment goat anti-mouse IgG, Fc_γ fragment specific secondary antibody labeled with the far-red fluorescent dye CF[®]640R.

Product Description

This product is prepared by labeling high quality F(ab')₂ fragment goat anti-mouse IgG, Fc_γ fragment specific, with the far-red fluorescent dye CF[®]640R. The antibody has been adsorbed against human, bovine and horse serum to minimize cross-reactivity. CF[®]640R is a better alternative to Cy[®]5, Alexa Fluor[®] 647 and other spectrally similar dyes. [Learn more about CF[®]640R.](#)

Advantages of CF[®]640R

- A novel rhodamine-based dye spectrally very similar to Cy[®]5 and Alexa Fluor[®] 647
- Much brighter than Cy[®]5 and at least as bright as Alexa Fluor[®] 647 when excited at 633 or 635 nm
- Exceptional photostability: far more photostable than Cy[®]5 and Alexa Fluor[®] 647
- More photostable than Atto 647N, a dye that has been extensively used in single molecule-based imaging
- Extremely water-soluble and pH-insensitive.

CF[®] Dyes offer exceptional brightness and photostability. For more information see our [CF[®] Dye technology page.](#)

Storage and Handling

Liquid format: Store at -20 °C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Liquid format antibodies contain 50% glycerol and will not freeze at -20 °C.

Lyophilized format: Store at -20 °C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Reconstitute antibodies in water using the indicated volumes below:

CF[®] Dye and biotin conjugates: add 0.5 mL dH₂O

HRP or DNP conjugates: add 1 mL dH₂O

Add the indicated volume of water directly to the vial containing the lyophilized antibody and mix gently to dissolve. Store reconstituted antibody at -20 °C and protect from light. Aliquot to avoid repeated freeze/thaw cycles. Alternatively, an equal volume of glycerol can be mixed with the reconstituted antibody so that it will remain liquid at -20 °C.

Optional: A preservative such as 0.05% sodium azide (final concentration) can be added to CF[®] Dye and biotin conjugates. Do not add sodium azide to HRP conjugates.

Note: Storage of the antibody for more than a day at final working dilution is not recommended.

CF is a registered trademark of Biotium, Inc.

Alexa Fluor is a registered trademark of Thermo Fisher Scientific. Cy Dye is a registered trademark of Cytiva.

References

Download a list of curated [CF[®] Dye references.](#)

Product attributes

Clonality	Polyclonal
Antibody type	Secondary
Antibody/conjugate formulation	Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution
Species reactivity	Mouse
Concentration	1 mg/mL (after reconstitution)
Host species	Goat
Antibody reactivity (target)	Mouse IgG, Fc
Cross adsorption	Bovine, Horse, Human
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.