

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, Fcγ 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, Fcγ 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](#).

This datasheet was generated on June 30, 2024 at 08:26:50 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf640r-goat-anti-mouse-igg-hl-fab2-fragment-fab2-fragment-fc%ce%b3-fragment-specific-highly-cross-adsorbed/>

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	Mouse IgG, Fc
Cross adsorption	Bovine, Horse, Human
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot