



Glowing products for science

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](#).

This datasheet was generated on December 24, 2025 at 09:50:26 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/>

Call us : [800-304-5357](tel:800-304-5357)

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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.