

GST Polyclonal Goat Antibody

Polyclonal goat anti-GST antibody labeled with our superior CF® Dyes.



Product attributes

Clonality	Polyclonal
Antibody type	Tag Antibody
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
Concentration	1 mg/mL
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Host species	Goat
Antibody reactivity (target)	GST

Product Description

This is a polyclonal goat anti-GST antibody labeled with our superior CF® Dyes. This antibody is designed to detect glutathione-S transferase (GST), a common expression tag used in molecular biology.

- Available in 5 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology

See our full selection of [anti-tag and anti-hapten antibody conjugates](#). View our full selection of [Secondary Antibodies](#), or search our catalog using our [Antibody Finder](#). Alternatively, you can view our [secondary antibody product listings](#) with catalog numbers. 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.

References

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 uL (100 ug)	20424	CF®488A Features
CF®594	593/615 nm	100 uL (100 ug)	20425	CF®594 Features
CF®640R	642/663 nm	100 uL (100 ug)	20426	CF®640R Features
CF®680	681/698 nm	50 uL (50 ug)	20981	CF®680 Features
CF®770	770/797 nm	50 uL (50 ug)	20982	CF®770 Features

This datasheet was generated on July 17, 2024 at 09:23:29 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/goat-anti-gst/>