

## 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
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.
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](#).

## GST Polyclonal Goat Antibody

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

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 dH<sub>2</sub>O

HRP or DNP conjugates: add 1 mL dH<sub>2</sub>O

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](#).

This datasheet was generated on February 14, 2025 at 04:16:42 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-gst/>