

Digoxigenin (DIG) Polyclonal Goat Antibody



Polyclonal goat anti-DIG antibody labeled with our superior CF® dyes and HRP.

Product Description

This is a polyclonal goat anti-DIG antibody labeled with our superior CF® dyes and HRP. This antibody recognizes the digoxigenin (DIG) hapten for detection of DIG conjugated proteins and nucleic acids.

- Available in 6 bright and photostable CF® dyes and HRP
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

See our full selection of [anti-tag and anti-hapten antibody conjugates](#).

Product attributes

Clonality	Polyclonal
Antibody type	Primary, Tag Antibody
Concentration	1 mg/mL (HRP conjugate), 2 mg/mL
Host species	Goat
Antibody reactivity (target)	DIG
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
Cross adsorption	Not cross-adsorbed
Secondary/tag antibody applications	ELISA, 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.

Goat Anti-DIG Tag

Conjugation	Ex/Em	Size	Catalog No.
CF®405S 250 uL (500 ug)	404/431 nm 20858-250uL	50 uL (100 ug)	20858-50uL
CF®488A 250 uL (500 ug)	490/515 nm 20859-250uL	50 uL (100 ug)	20859-50uL
CF®568 250 uL (500 ug)	562/583 nm 20860-250uL	50 uL (100 ug)	20860-50uL
CF®594 250 uL (500 ug)	593/614 nm 20861-250uL	50 uL (100 ug)	20861-50uL
CF®640R 250 uL (500 ug)	642/662 nm 20862-250uL	50 uL (100 ug)	20862-50uL
CF®680R 250 uL (500 ug)	680/701 nm 20863-250uL	50 uL (100 ug)	20863-50uL
HRP	N/A	100 uL (100 ug)	20864-100uL

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](#).

This datasheet was generated on June 3, 2026 at 03:05:35 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf-dye-hrp-dig-tag-polyclonal-goat-antibody/>