

Digoxigenin (DIG) Polyclonal Goat Antibody

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



Product attributes

Clonality	Polyclonal
Antibody type	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 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](#). 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.
CF®405S	404/431 nm	50 uL (100 ug)	20858-50uL
		250 uL (500 ug)	20858-250uL
CF®488A	490/515 nm	50 uL (100 ug)	20859-50uL
		250 uL (500 ug)	20859-250uL
CF®568	562/583 nm	50 uL (100 ug)	20860-50uL
		250 uL (500 ug)	20860-250uL
CF®594	593/614 nm	50 uL (100 ug)	20861-50uL
		250 uL (500 ug)	20861-250uL
CF®640R	642/662 nm	50 uL (100 ug)	20862-50uL
		250 uL (500 ug)	20862-250uL
CF®680R	680/701 nm	50 uL (100 ug)	20863-50uL
		250 uL (500 ug)	20863-250uL
HRP	N/A	100 uL (100 ug)	20864-100uL