

# Digoxigenin polyclonal antibody

Species independent all-purpose immuno-tag.

Digoxigenin (DIG) is a steroid found exclusively in the flowers and leaves of the plants *Digitalis purpurea* and *Digitalis lanata*. DIG is the steroid component of digoxin, a cardiac glycoside, and can be conjugated onto a variety of biomolecules (proteins, nucleic acids) to be detected in further assays (e.g. ELISA). These antibodies can be labeled with dyes, enzymes or fluorescence, directly or secondarily, for visualization and detection.

DIG has an advantage over biotin due to the fact that biotin is an endogenous molecule of living cells and plays a key role in many reactions, which can sometimes lead to false positive results. Because the blossoms and leaves of *Digitalis* plants are the only natural sources of digoxigenin, binding of the anti-digoxigenin antibody in other biological material is not a consideration.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-ABS302-0001	1mg
-----------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	ELISA, ISH (in situ hybridization), TUNEL, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Sheep
Immunogen	Digoxin
Purity Detail	Antigen affinity purified.
Source	Purified from sheep serum.
Species Reactivity	Species independent
Specificity	Recognizes Digoxin and Digoxigenin.
Technical Info / Product Notes	This product is unconjugated but if you desire a specific conjugation please contact technical support for lead time and pricing. No minimum quantity required.
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)