

# Green fluorescent protein polyclonal antibody

Green Fluorescent Protein (GFP) is responsible for bioluminescence in the jellyfish *Aequorea victoria*. This protein has quickly become a powerful research tool for assessing gene expression and subcellular protein distribution in fixed or living cells. GFP is excited by and brightly fluoresces when exposed to UV or blue light. This feature makes it ideal as a marker for use in fluorescence microscopy, cytometry, tagging fusion proteins, and assaying transcriptional regulation from gene promoters in vivo. Numerous GFP variants with enhanced and shifted emission spectra (blue, green, and yellow) have been developed through aa substitutions at specific residues.

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

Citations: 6

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-210-1990	100µl
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** GFP

**Application** ICC, WB

**Application Notes** Detects a band of ~28 kDa by Western blot.

**Formulation** Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 26-39 (H26KFSVSGEGEGDAT39) of the jellyfish Aequorea victoria GFP (green fluorescent protein). This sequence is completely conserved between GFP and the modified and enhanced Blue, Green and Yellow FP variants.

**Purity Detail** Epitope-affinity purified.

**Recommendation Dilutions/Conditions** Immunocytochemistry (1:200)Western Blot (0.5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Species independent

**Specificity** Recognizes GFP and GFP fusion proteins.

**UniProt ID** P42212

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

Last modified: May 29, 2024



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)