

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-210-199-R100	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](#).



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)