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 as 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 NotesDetects 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



info-

eu@enzolifesciences.com