

# Green Fluorescent Protein monoclonal antibody (1E4)

The green fluorescent protein (GFP) is derived from the jellyfish *Aequorea victoria*, that fluoresces green when exposed to blue light. GFP is a frequently used reporter for gene expression in recombinant DNA and cell transfection systems.

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

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-SAB-500-E	100µg
---------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name GFP

Application ICC, IHC, IP, WB

Application Notes Detects a band of ~27kDa by Western blot.

Clone 1E4

Formulation Liquid. In PBS containing 50% glycerol.

Gene/Protein Identifier JQ1514 (PIR)

Host Mouse

Immunogen Recombinant GFP protein (Aequorea victoria).

Isotype IgG2b

Purity Detail Protein A affinity purified.

**Recommendation** Immunocytochemistry (2µg/ml)Immunohistochemistry (10µg/ml)Immunoprecipitation  
**Dilutions/Conditions** (5µg/ml)Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Species independent

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)