

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 Dilutions/Conditions	Immunocytochemistry (2µg/ml)Immunohistochemistry (10µg/ml)Immunoprecipitation (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)