

AbX™

Green Fluorescent Protein (GFP) Polyclonal Antibody

Catalog Number A013-25UG



ARBOR
ASSAYS

FEATURES

- Rabbit Antibody to wild type Green Fluorescent Protein (GFP)
- Affinity Purified on a GFP column
- Reacts with all GFP variants

INTRODUCTION

Green fluorescent protein (GFP), a 27 kDa protein derived from the jellyfish *Aequorea victoria*, emits green light (emission peak 509 nm) when excited by blue light (excitation peak 395 nm). GFP has become an invaluable tool in cell biology research, since its intrinsic fluorescence can be visualized in living cells. GFP fluorescence is stable under fixation conditions and suitable for a variety of applications. GFP has been widely used as a reporter for gene expression, enabling researchers to visualize and localize GFP-tagged proteins within living cells without the need for chemical staining. Other applications of GFP include assessment of protein-protein interactions through the yeast two hybrid system and measurement of distance between proteins through fluorescence energy transfer (FRET) protocols. GFP is used to measure single cell metastasis and successful proliferation of stem cells.

FORM:	100 mM Sodium Phosphate, 150 mM Sodium Chloride, 0.09% Na Azide, pH 7.2.	
PURITY:	Affinity purified on a GFP column. Single IgG peak	
CONCENTRATION:	100 µg/mL	
STORAGE:	4°C for short term, up to 24 months at -20°C.	
IMMUNOGEN:	Highly purified native GFP from <i>Aequorea victoria</i> .	
SPECIFICITY:	It is reactive with GFP and GFP fusion proteins from native and recombinant sources. Recognizes all variants of GFP.	
USES:	For Immunoblotting	Dilute 1:1,000 or greater
COUNTRY OF ORIGIN:	USA	

FOR RESEARCH USE ONLY