

PE/XFD 700 Goat Anti-Mouse IgG (H+L) Antibody

Catalog Number: 50323

Unit Size: 200 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	~0.2 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Mouse
Class	Secondary
Clonality	Polyclonal
Host	Goat

Biological Properties

Stabilizer	0.09% sodium azide, 0.2% (w/v) BSA
Appearance	Liquid
Preparation	Goat anti-mouse IgG (H+L) is produced in goat with pooled total mouse IgG. The antibody is conjugated with PE/XFD 700 under optimal conditions.
Application	Flow Cytometry (FACS), IF, IHC
Recommended Dilutions	Suggested dilutions are only guidelines; users should titrate the product for their specific assay using appropriate controls

Application	Recommended dilution
Flow Cytometry (FACS)	1-5 µg/mL
IF	2 µg/mL
IHC	1-10 µg/mL

Spectral Properties

Conjugate	PE (phycoerythrin)-Alexa Fluor 700
-----------	------------------------------------

Excitation Wavelength 565 nm

Emission Wavelength 721 nm

Applications

PE/XFD 700 Goat Anti-Mouse IgG (H+L) Antibody is a high-performance secondary antibody designed to bind both heavy and light chains of mouse immunoglobulin G (IgG). Conjugated to PE/XFD 700, this antibody delivers exceptionally bright far red fluorescence with excitation/emission maxima of ~565/721 nm. The PE/XFD 700 fluorophore is prepared using AAT Bioquest's proprietary labeling technology to ensure superior signal intensity compared to similar antibodies from other sources. This antibody is ideal for flow cytometry, fluorescence microscopy, ELISA, and other immunoassays requiring sensitive and specific detection of mouse IgG. Its compatibility with 561 nm yellow-green laser lines and common bandpass filters (e.g., 730/45 nm) ensures seamless integration into standard fluorescence detection workflows.