

Purified Goat Anti-human IgG Fc (H+L) Antibody

Catalog Number: 50273

Unit Size: 1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Clonality	Monoclonal
Host	Goat
Immunogen	igg

Chemical Properties

Molecular Weight	~150000
------------------	---------

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Soluble In	PBS
Application	Flow Cytometry (FACS)

Applications

AAT Bioquest's anti-human secondary antibodies have well-characterized specificity for human immunoglobulins and are useful in the detection, sorting or purification of its specified target. They can also provide greater sensitivity through signal amplification as multiple secondary antibodies can bind to a single primary antibody. This antibody was purified through affinity chromatography.