

**PE-Cy5.7 Mouse Anti-human IgG (Fab)
Antibody *4A11, monoclonal***Catalog number: V1031485
Unit size: 100 tests**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), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Secondary
Clonality	Monoclonal
Host	Mouse
Immunogen	IgG (Fab)
Clone	4A11
Conjugate	PE-Cy5.7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE-Cy5.7 under optimal conditions
Application	FC (QC TESTED)

Applications

Detects the antigen-binding fragment of Immunoglobulin G.