

FITC Mouse Anti-human IgG (Fc) Antibody
EM-07, monoclonalCatalog number: V1031500
Unit size: 0.1 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), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Secondary
Clonality	Monoclonal
Host	Mouse
Immunogen	IgG (Fc)
Clone	EM-07
Conjugate	FITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions
Application	FC (QC TESTED)

Applications

High affinity immunoglobulin γ Fc receptor I (IgG Fc receptor I) is a 49 kDa transmembrane protein that can be found in the clathrin-coated endocytic vesicle membrane, plasma membrane and early endosome membrane of cells. It is alternatively called Fc- γ RIA (IgG Fc receptor I) and Fc- γ RI (IgG Fc receptor I). In humans, Fc- γ RIA plays a role in the upregulation of protein tyrosine kinase activity. It is the subject of comprehensive study due to the fact that it plays a role in the Fc- γ receptor signaling pathway involved in phagocytosis and interferon- γ -mediated signaling pathway. Sequencing of Fc- γ RIA has exemplified it contains 5 conserved structural units: extracellular, cytoplasmic, Ig-like C2-type 1, Ig-like C2-type 2 and Ig-like C2-type 3 domain. It is thought to be essential to immune response. It binds to IgG. And it plays an important role in organismal processes, in particular, immune response, antigen processing and presentation of exogenous peptide antigen via MHC class I, TAP-dependent and receptor-mediated endocytosis.