

Purified Mouse Anti-human PAG1 Antibody
MEM-255, monoclonal, Cross AdsorbedCatalog number: V1031965
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	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	PAG1
Clone	MEM-255

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse, rat, cow serum and then conjugated with under optimal conditions
Application	FC, WB, IHC(P)

Applications

Phosphoprotein associated with glycosphingolipid-enriched microdomains 1 (alternatively called Transmembrane adapter protein PAG, Transmembrane phosphoprotein Cbp or Csk-binding protein) is a transmembrane protein with a molecular weight of 47 kDa, found in the membrane raft, integral component of membrane and plasma membrane of cells. Sequencing of transmembrane adapter protein PAG has shown evidence that it contains 2 types of conserved structural units: extracellular and cytoplasmic domain. Transmembrane adapter protein PAG aids in T cell activation. It is the subject of intensive research stemming from the fact that it is involved with the T cell receptor signaling pathway. Transmembrane adapter protein PAG is thought to be essential to organismal processes, for example, intracellular signal transduction, adaptive immune response and signal transduction. It acts to negatively regulate T cell activation. Transmembrane adapter protein PAG binds with SH2 domain and S-palmitoyl cysteine.