

PE Mouse Anti-human ARHGEF4 Antibody
ARHGEF-08, monoclonalCatalog number: V103135
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	ARHGEF4
Clone	ARHGEF-08
Conjugate	PE

Biological Properties

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

Applications

p guanine nucleotide exchange factor 4 is a protein with a molecular weight of 79 kDa, located in the ruffle membrane and cytosol of cells. In humans, p guanine nucleotide exchange factor 4 plays an important role in small GTPase mediated signal transduction. It has been associated with key functions like Rac guanyl-nucleotide exchange factor and guanyl-nucleotide exchange factor activity. Sequencing of p guanine nucleotide exchange factor 4 has exemplified it contains 3 types of conserved structural units: SH3, DH and PH domain. p guanine nucleotide exchange factor 4 enhances apoptotic process, and furthermore, binds with protein domain specific. p guanine nucleotide exchange factor 4 is the subject of comprehensive study because of the fact that it is a member of the G protein-coupled receptor signaling pathway. It plays an important role in organismal processes, for example, intracellular signal transduction, lamellipodium assembly and filopodium assembly.