

**Purified Mouse Anti-human PDE8A
Antibody *EM-52, monoclonal***

Catalog number: V1031995

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	PDE8A
Clone	EM-52

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB

Applications

High affinity cAMP-specific and IBMX-insensitive 3',5'-cyclic phosphodiesterase 8A is a protein with a molecular weight of 93 kDa, located in the extracellular exosome and cytosol of cells. In humans, High affinity cAMP-specific and IBMX-insensitive 3',5'-cyclic phosphodiesterase 8A is the subject of comprehensive research in part because it is involved with the G protein-coupled receptor signaling pathway. It has been closely linked to critical functions like 3',5'-cyclic-AMP phosphodiesterase and 3',5'-cyclic-GMP phosphodiesterase activity. Sequencing of high affinity cAMP-specific and IBMX-insensitive 3',5'-cyclic phosphodiesterase 8A has exemplified it contains 3 types of conserved structural units: PAC, PAS and PDEase domain. High affinity cAMP-specific and IBMX-insensitive 3',5'-cyclic phosphodiesterase 8A has been found to be involved in transcription, DNA-templated transcription. It is an integral part of organismal processes, for example, cellular response to epidermal growth factor stimulus and cAMP catabolic process. High affinity cAMP-specific and IBMX-insensitive 3',5'-cyclic phosphodiesterase 8A is a positive regulator of ERK1 and ERK2 cascade and protein phosphorylation. It also plays a role in the downregulation of hydrogen peroxide-induced cell death and cell death. It binds with metal ion and kinase.