

**Purified Mouse Anti-human PRKAR2A
Antibody *Hs-36, monoclonal***

Catalog number: V1032035

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	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	PRKAR2A
Clone	Hs-36

Biological Properties

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

Applications

cAMP-dependent protein kinase type II- α regulatory subunit is a 46 kDa protein that can be located in the nucleotide-activated protein kinase complex, cytosol and centrosome of cells. In humans, cAMP-dependent protein kinase type II- α regulatory subunit is an inhibitor of cAMP-dependent protein kinase activity. cAMP-dependent protein kinase type II- α regulatory subunit has been closely linked to critical functions such as cAMP-dependent protein kinase regulator and cAMP-dependent protein kinase inhibitor activity. It reacts with protein kinase A catalytic subunit, cAMP 1 and cAMP. cAMP-dependent protein kinase type II- α regulatory subunit is essential to protein kinase A signaling and cAMP-mediated signaling. It aids in organismal processes, for instance, renal water homeostasis, intracellular signal transduction and activation of protein kinase A activity.