## PIP<sub>2</sub> monoclonal antibody (KT10)

Metabolism of inositol phospholipids by intracellular signaling mediators is fundamental to signal transduction in eukaryotic cells. PI-4,5-P2 (PIP2) can be synthesized by phosphorylation of PI-4-P by type I phosphatidylinositol phosphate kinase (PIP5K I), or phosphorylation of PI-5-P by type II PIPK (PIP4K II). PI-4,5-P2 regulation of cellular calcium levels involves its hydrolysis by Phospholipase C (PLC) to produce inositol 1,4,5-triphosphate (IP3) and diacylglycerol (DAG), which serve as second messengers in the import of calcium via IP3-sensitive ion channels and in the activation of PKC, respectively.

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

**Ordering Information** 

**Order Online** »

ADI-915-052-020

20µl

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name Phosphatidylinositol 4,5-bisphosphate, PtdIns(4,5)P2

Application ELISA, ICC

Clone KT10

Formulation Liquid. Ascites.

**Host** Mouse

**Immunogen** Native PIP2 from bovine spinal cord.

lsotype lgG2b

**Source** Purified from mouse ascites.

Species Reactivity Species independent

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

