

# 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-P<sub>2</sub> (PIP<sub>2</sub>) 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-P<sub>2</sub> regulation of cellular calcium levels involves its hydrolysis by Phospholipase C (PLC) to produce inositol 1,4,5-triphosphate (IP<sub>3</sub>) and diacylglycerol (DAG), which serve as second messengers in the import of calcium via IP<sub>3</sub>-sensitive ion channels and in the activation of PKC, respectively.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 3

[View Online »](#)

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.
Isotype	IgG2b
Source	Purified from mouse ascites.
Species Reactivity	Species independent
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)