

PIP₂ 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.

Isotype IgG2b

Source Purified from mouse ascites.

Species Reactivity Species independent

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



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)