

PRODUCT INFORMATION



PtdIns-(3,4,5)-P₃ Binding Protein Item No. 10009817

Overview and Properties

Synonyms:	Cyth3, Cytohesin 3, GRP1, GRP1 PH Domain, PIP ₃ , PIP ₃ [3',4',5'] Binding Protein, PI(3,4,5)-P ₃ , PI(3,4,5)P ₃ Binding Protein, Phosphatidylinositol-3,4,5-triphosphate
Source:	Recombinant human N-terminal GST-tagged PIP ₃ expressed in <i>E. coli</i>
Amino Acids:	264-381 (PH domain)
Uniprot No.:	O43739
Molecular Weight:	~41.7 kDa (15.4 kDa native; 26.3 kDa as provided)
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥95%
Supplied in:	A solution in PBS, pH 7.2, containing 20% glycerol
Protein Concentration:	1.0 mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images

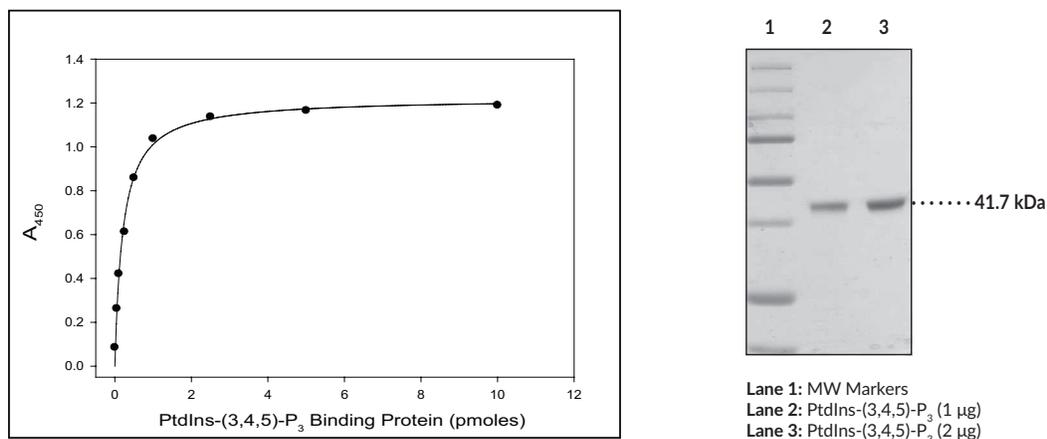


Figure 1: PtdIns-(3,4,5)-P₃-biotinamide (10 pmol/well) was applied to a streptavidin-coated, 96-well plate and was detected using PtdIns-(3,4,5)-P₃ binding protein. Colorimetric readout at 450 nm was accomplished using a HRP-conjugated anti-GST antibody.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/19/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Pleckstrin homology (PH) domains all contain a seven-standard β -sandwich that allows binding to various species of phosphatidylinositol (PtdIns) phosphates.¹ Most PH domains bind to PtdIns phosphates with weak affinity and low specificity. However, a small subclass binds specifically and with high-affinity for certain PtdIns phosphates. PtdIns-(3,4,5)-P₃ Binding Protein contains a highly specific PH domain that recognizes and binds PtdIns-(3,4,5)-P₃ (see Figure 1).² PtdIns phosphates represent a small percentage of total membrane phospholipids. However, they play a critical role in the generation and transmission of cellular signals.^{3,4} PtdIns-(4,5)-P₂ can be phosphorylated by phosphoinositide (PI)-3-kinase to make PtdIns-(3,4,5)-P₃ which initiates an intricate signalling cascade that has been implicated in cancer.⁵ PtdIns-(3,4,5)-P₃ binding protein can be used in *in vitro* assays for the detection of PtdIns-(3,4,5)-P₃ in PI₃-kinase and PTEN phosphatase assays. Cayman's PtdIns-(3,4,5)-P₃ Binding Protein is isoform 2, which differs from isoform 1 by the absence of a glycine residue at position 277.

References

1. Lemmon, M.A. and Ferguson, K.M. *Biochem. Soc. Trans.* **29**(4), 377-384 (2001).
2. Lemmon, M.A. *Traffic* **4**, 201-213 (2003).
3. Exton, J.H. *Annu. Rev. Pharmacol. Toxicol.* **36**, 481-509 (1996).
4. Majerus, P.W. *Annu. Rev. Biochem.* **61**, 225-250 (1992).
5. Vivanco, I. and Sawyers, C.L. *Nature Reviews Cancer* **2**, 489-501 (2002).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM