

PRODUCT INFORMATION



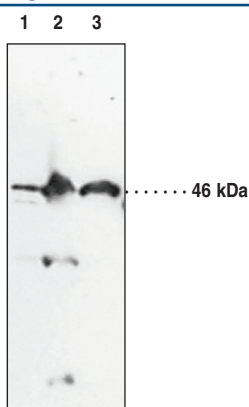
NAPE-PLD (N-Term) Polyclonal Antibody

Item No. 10306

Overview and Properties

Contents:	This vial contains 500 µl of peptide-affinity purified antibody.
Synonyms:	N-Acyl-Phosphatidylethanolamine-Hydrolysing Phospholipase D
Immunogen:	Peptide from the N-terminal region of human NAPE-PLD
Species Reactivity:	Human, mouse, rat, and bovine NAPE-PLD (N-Term)
Uniprot No.:	Q61Q20
Form:	Liquid
Storage:	-20°C (as supplied)
Storage Buffer:	TBS, pH 7.4, 50% glycerol, 0.1% BSA, and 0.02% sodium azide
Stability:	≥1 year
Host:	Rabbit
Application:	Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image(s)



Lane 1: Human Cerebellum Supernatant (30 µg)
Lane 2: Mouse Brain Homogenate (50 µg)
Lane 3: Mouse Brain High-Density Membrane (30 µg)

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, 03/04/2016

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

N-acylethanolamines (NAEs) are involved in diverse biological processes such as inflammatory regulation, apoptosis, and tissue degeneration.¹ In animals, NAEs are mainly biosynthesized *via* a membrane phospholipid-dependent pathway, which is the enzymatic hydrolysis of N-acyl-phosphatidylethanolamine (NAPE). The enzyme catalyzing this reaction is a phospholipase D subtype selective for NAPE named N-acyl-phosphatidylethanolamine-hydrolysing phospholipase D (NAPE-PLD). It has been cloned from mouse, rat, and human and is 393-396 amino acids in length, with an estimated molecular weight of 46 kDa. Both NAPE-PLD mRNA and protein activity have been detected in a wide range of tissues with the highest levels in brain, kidney, and testis.² In rat, NAPE-PLD activity in the brain is low in neonates and is 15-fold higher in adults, whereas the activity remains constant in the heart during development.³

References

1. Hansen, H.S., Moesgaard, B., Petersen, G., *et al.* Putative neuroprotective actions of N-acyl-ethanolamines. *Pharmacology & Therapeutics* **95**, 119-126 (2002).
2. Okamoto, Y., Morishita, J., Tsuboi, K., *et al.* Molecular characterization of a phospholipase D generating anandamide and its congeners. *J. Biol. Chem.* **279**(7), 5298-5305 (2004).
3. Moesgaard, B., Petersen, G., Jaroszewski, J.W., *et al.* Age dependent accumulation of N-acyl-ethanolamine phospholipids in ischemic rat brain: A ³¹P NMR and enzyme activity study. *J. Lipid Res.* **41**, 985-990 (2000).

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