

PRODUCT INFORMATION



ABHD4 (human, recombinant)

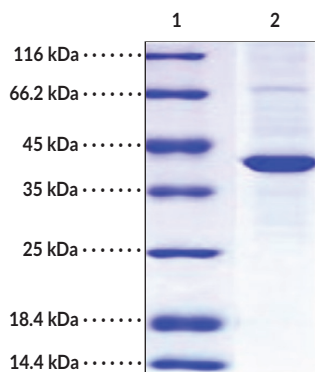
Item No. 44970

Overview and Properties

Synonyms: Abhydrolase Domain-containing Protein 4, α/β -Hydrolase 4, α/β Hydrolase Domain-containing Protein 4, (Lyso)-N-Acylphosphatidylethanolamine Lipase
Source: Recombinant human N-terminal His-tagged ABHD4 expressed in insect cells
Amino Acids: 1-342 (full length)
Uniprot No.: Q8TB40
Molecular Weight: 41 kDa
Storage: -80°C (as supplied)
Stability: ≥ 1 year
Purity: $\geq 82\%$ estimated by SDS-PAGE
Supplied in: Lyophilized from sterile 50 mM Tris, with 100 mM sodium chloride, 10% glycerol, and pH 8.0
Endotoxin Testing: < 1.0 EU/ μ g, determined by the LAL endotoxin assay
Protein Concentration: *batch specific* mg/ml

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

Image



Lane 1: MW Markers
Lane 2: ABHD4

SDS-PAGE Analysis of ABHD4. This protein has a calculated molecular weight of 41 kDa.

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, 12/15/2025

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

α/β -Hydrolase domain-containing protein 4 (ABHD4) is a serine hydrolase and lysophospholipase/phospholipase B that hydrolyzes N-acyl phosphatidylethanolamines (NAPEs) and lyso-NAPEs.^{1,2} It is expressed in the brain, spinal cord, testes, liver, kidney, and heart.² *Abhd4*^{-/-} mice exhibit decreased brain levels of various N-acyl phospholipids, including plasmalogen-based lyso-NAPEs, N-acyl lysophosphatidylserines, and glycerophospho-N-acylethanolamines (GP-NAEs).¹ Cayman's ABHD4 (human, recombinant) protein consists of 360 amino acids and has a calculated molecular weight of 41 kDa.

References

1. Lee, H.-C., Simon, G.M., and Cravatt, B.F. ABHD4 regulates multiple classes of N-acyl phospholipids in the mammalian central nervous system. *Biochemistry* **54(15)**, 2539-2549 (2015).
2. Simon, G.M. and Cravatt, B.F. Endocannabinoid biosynthesis proceeding through glycerophospho-N-acyl ethanolamine and a role for α/β -hydrolase 4 in this pathway. *J. Biol. Chem.* **281(36)**, 26465-26472 (2006).

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