

PRODUCT INFORMATION



SMPDL3A (human, recombinant; His-tagged)

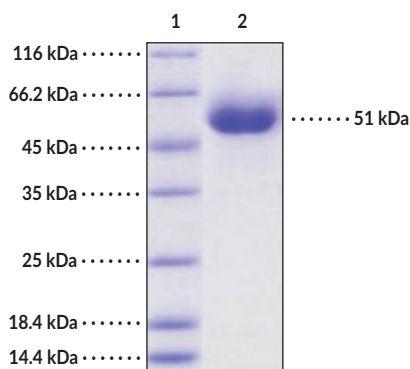
Item No. 40468

Overview and Properties

Synonyms: Acid Sphingomyelinase-like Phosphodiesterase 3A, ASML3A, ASM-like Phosphodiesterase 3A, Sphingomyelin Phosphodiesterase Acid-like 3A
Source: Recombinant C-terminal His-tagged SMPDL3A expressed in HEK293 cells
Amino Acids: 1-453
Molecular Weight: 50.3 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥95% estimated by SDS-PAGE
Supplied in: Lyophilized from sterile PBS, pH 7.4
Endotoxin Testing: <1.0 EU/μg, determined by the LAL endotoxin assay

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

Image



Lane 1: MW Markers
Lane 2: SMPDL3A

SDS-PAGE Analysis of SMPDL3A. This protein has a calculated molecular weight of 50.3 kDa. It has an apparent molecular weight of approximately 51 kDa by SDS-PAGE under reducing conditions due to glycosylation.

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, 02/01/2024

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

Sphingomyelin phosphodiesterase acid-like 3A (SMPDL3A) is a zinc-dependent nucleotide phosphoesterase and member of the acid sphingomyelinase-like enzyme family.¹ It is composed of an N-terminal catalytic domain and a C-terminal domain, which contains an asparagine located at position 356 (Asn356). The N-glycosylation of Asn356 is required for SMPDL3A stability, enzymatic activity, and cellular secretion.^{1,2} SMPDL3A is expressed in, and secreted from, osteoclasts, adipocytes, astrocytes, and macrophages and reduces local inflammation induced by nucleotide tri- or diphosphates.^{1,3} Its intracellular levels and secretion are increased by cAMP, cholesterol, and liver X receptor (LXR) activation, and it localizes to the lysosome, which provides the acidic environment required for its optimal catalytic activity.^{1,3} SMPDL3A knockout decreases cell proliferation, colony formation, and migration, as well as induces cell cycle arrest in the G₀/G₁ phase and apoptosis in HepG2 hepatocellular carcinoma cells.⁴ *In vivo*, SMPDL3A knockout in HepG2 cells reduces tumor volume and weight in a HepG2 xenograft mouse model. SMPDL3A levels are increased in patients with hepatocellular carcinoma (HCC) and negatively correlated with prognosis. Cayman's SMPDL3A (human, recombinant; His-tagged) protein consists of 442 amino acids and has a calculated molecular weight of 50.3 kDa. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 51 kDa due to glycosylation.

References

1. Gorelik, A., Illes, K., Superti-Furga, G., *et al.* Structural basis for nucleotide hydrolysis by the acid sphingomyelinase-like phosphodiesterase SMPDL3A. *J. Biol. Chem.* **291**(12), 6376-6385 (2016).
2. Traini, M., Kumaran, R., Thaysen-Andersen, M., *et al.* N-glycosylation of human sphingomyelin phosphodiesterase acid-like 3A (SMPDL3A) is essential for stability, secretion and activity. *Biochem. J.* **474**(7), 1071-1092 (2017).
3. Traini, M., Quinn, C.M., Sandoval, C., *et al.* Sphingomyelin phosphodiesterase acid-like 3A (SMPDL3A) is a novel nucleotide phosphodiesterase regulated by cholesterol in human macrophages. *J. Biol.Chem.* **289**(47), 32895-32913 (2014).
4. Zhang, Y., Chen, W., Cheng, X., *et al.* Sphingomyelin phodiesterase acid-like 3A promotes hepatocellular carcinoma growth through the enhancer of rudimentary homolog. *Front. Oncol.* **12**, 852765 (2022).

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