

PRODUCT INFORMATION



ABHD14B (human, recombinant)

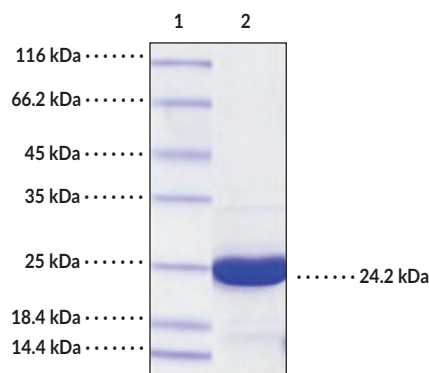
Item No. 44971

Overview and Properties

| | |
|--------------------------|---|
| Synonyms: | Abhydrolase Domain-containing 14B, α/β -Hydrolase Domain-containing Protein 14B, CCG1-interacting Factor B, CCG1/TAFII250-interacting Factor B, CCG1/TAF(II)250-interacting Factor B, CIB, Putative Protein-lysine Deacylase ABHD14B |
| Source: | Recombinant human N-terminal His-tagged ABHD14B expressed in <i>E. coli</i> |
| Amino Acids: | 1-210 |
| Uniprot No.: | Q96IU4 |
| Molecular Weight: | 24.2 kDa |
| Storage: | -80°C (as supplied) |
| Stability: | ≥ 1 year |
| Purity: | $\geq 96\%$ estimated by SDS-PAGE |
| Supplied in: | Supplied as sterile PBS, pH 7.4, with 20% glycerol, 50 mM arginine, and 0.1% polysorbate 20 Protein |
| Concentration: | <i>batch specific</i> |

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

Image



Lane 1: MW Markers
Lane 2: ABHD14B

SDS-PAGE Analysis of ABHD14B. ABHD14B has a calculated molecular weight of 24.2 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, 12/12/2025

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 14B (ABHD14B) is a lysine deacetylase and member of the metabolic serine hydrolase family.^{1,2} It is composed of an ABHD fold that contains the Ser-His-Asp catalytic triad. ABHD14B is expressed in kidney, liver, and lung and localizes to the cytosol and nucleus. It transfers an acetyl group from acetylated lysines on proteins, such as histones, to CoA to produce acetyl-CoA.^{1,3} ABHD14B is involved in the regulation of glycolysis and the citric acid cycle, as well as viral transcription from hepatitis B virus (HBV) covalently closed circular DNA (cccDNA).^{2,3} Decreased expression of ABHD14B is associated with poor prognosis in patients with breast cancer.⁴ Cayman's ABHD14B (human, recombinant) protein consists of 225 amino acids and has a calculated molecular weight of 24.2 kDa. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 24 kDa.

References

1. Rajendran, A., Vaidya, K., Mendoza, J., *et al.* Functional annotation of ABHD14B, an orphan serine hydrolase enzyme. *Biochemistry* **59**(2), 183-196 (2020).
2. Rajendran, A., Soory, A., Khandelwal, N., *et al.* A multi-omics analysis reveals that the lysine deacetylase ABHD14B influences glucose metabolism in mammals. *J. Biol. Chem.* **298**(7), 102128 (2022).
3. Sekiba, K., Miyake, N., Miyakawa, Y., *et al.* CRISPR-mediated proximity labeling unveils ABHD14B as a host factor to regulate HBV cccDNA transcriptional activity. *Hepatology* **9**(8), e0757 (2025).
4. Aushev, V.N., Gopalakrishnan, K., Teitelbaum, S.L., *et al.* Tumor expression of environmental chemical-responsive genes and breast cancer mortality. *Endocr. Relat. Cancer* **26**(12), 843-851 (2019).

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