

PRODUCT INFORMATION



PDE3A (human, recombinant)

Item No. 32041

Overview and Properties

Synonyms: CGI-PDE A, cGMP-inhibited 3',5'-Cyclic Phosphodiesterase A, PDE3A, Phosphodiesterase 3A, CGMP-inhibited

Source: Recombinant human N-terminal His-GST-tagged PDE3A expressed in insect cells

Amino Acids: 669-1141

Uniprot No.: Q14432

Molecular Weight: 81.7 kDa

Storage: -80°C (as supplied)

Stability: ≥1 year

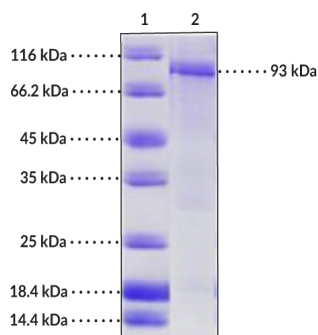
Purity: ≥90% estimated by SDS-PAGE

Supplied in: Lyophilized from sterile 20 mM Tris, pH 7.4, with 500 mM sodium chloride and 10% glycol, 3 mM DTT

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: PDE3A

SDS-PAGE Analysis of PDE3A. This protein consists of 710 amino acids and has a calculated molecular weight of 81.7 kDa. By SDS-PAGE, under reducing conditions, the molecular mass of the protein is 93 kDa due to apparent post-translational modifications.

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, 06/10/2020

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

Phosphodiesterase 3A (PDE3A) is a cyclic nucleotide phosphodiesterase that hydrolyzes cGMP and cAMP.^{1,2} It is a 125 kDa protein composed of an N-terminal domain important for localization, a linker region, and a conserved C-terminus that contains the catalytic domain.² Alternative splicing produces three isomers of PDE3A, PDE3A1, PDE3A2, and PDE3A3, that vary in N-terminal hydrophobic membrane association region (NHR) composition but all contain the same conserved C-terminal catalytic domain. PDE3A is found in the heart, vascular smooth muscle, platelets, and oocytes.³ Gain-of-function mutations in *PDE3A* have been found in individuals with Mendelian hypertension and brachydactyly type E.⁴ *PDE3A* expression is decreased in drug-resistant non-small cell lung cancer (NSCLC) cells, and overexpression of *PDE3A* restores sensitivity to cisplatin (Item No. 13119) in cisplatin-resistant A549 cells.⁵ Female *Pde3a* knockout mice exhibit inhibited oocyte maturation and complete sterility.⁶ *Pde3a* knockout mice also exhibit increased survival compared with *Pde3a* wild-type mice in a model of collagen and epinephrine-induced pulmonary thromboembolism. Cayman's PDE3A (human, recombinant) protein consists of 710 amino acids and has a calculated molecular weight of 81.7 kDa. By SDS-PAGE, under reducing conditions, the molecular mass of the protein is 93 kDa due to apparent post-translational modifications.

References

1. Kütthe, A., Eckel, H., Stief, C.G., *et al.* Molecular biological characterization of phosphodiesterase 3A from human corpus cavernosum. *Chem. Biol. Interact.* **119-120**, 593-598 (1999).
2. Vandeput, F., Szabo-Fresnais, N., Ahmad, F., *et al.* Selective regulation of cyclic nucleotide phosphodiesterase PDE3A isoforms. *Proc. Natl. Acad. Sci. USA* **110(49)**, 19778-19783 (2013).
3. Lugnier, C. Cyclic nucleotide phosphodiesterase (PDE) superfamily: A new target for the development of specific therapeutic agents. *Pharmacol. Ther.* **109(3)**, 366-398 (2006).
4. Maass, P.G., Aydin, A., Luft, F.C., *et al.* *PDE3A* mutations cause autosomal dominant hypertension with brachydactyly. *Nat. Genet.* **47(6)**, 647-653 (2015).
5. Tian, F.-M., Zhong, C.-Y., Wang, X.-N., *et al.* PDE3A is hypermethylated in cisplatin resistant non-small cell lung cancer cells and is a modulator of chemotherapy response. *Eur. Rev. Med. Pharmacol. Sci.* **21(11)**, 2635-2641 (2017).
6. Begum, N., Shen, W., and Manganiello, V. Role of PDE3A in regulation of cell cycle progression in mouse vascular smooth muscle cells and oocytes: Implications in cardiovascular diseases and infertility. *Curr. Opin. Pharmacol.* **11(6)**, 725-729 (2011).

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