

PRODUCT INFORMATION



CdnP (*Mycobacterium tuberculosis* strain ATCC 25618/H37Rv recombinant)

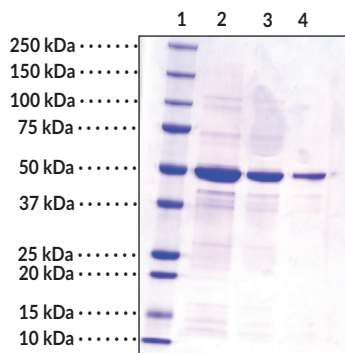
Item No. 22809

Overview and Properties

Synonyms:	Bifunctional Oligoribonuclease and PAP Phosphatase NrnA, 3'(2'),5'-Bisphosphate Nucleotidase, CnpB, Cyclic di-NMP Phosphodiesterase, PAP Phosphatase, 3'-Phosphoadenosine 5'-phosphate Phosphatase, Rv2837c
Source:	Active recombinant <i>M. tuberculosis</i> strain ATCC 25618/H37Rv N-terminal Trx-, His-, and S-tagged CdnP expressed in <i>E. coli</i>
Amino acids:	1-336 (full length)
Uniprot No.:	P71615
Molecular Weight:	52.71 kDa
Storage:	-80°C (as supplied); avoid freeze/thaw cycles by storing protein in aliquots
Stability:	≥1 year
Purity:	batch specific (≥70% estimated by SDS-PAGE)
Supplied in:	50 mM HEPES, pH 8.0, with 150 mM sodium chloride
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: CdnP (5 µg)
Lane 3: CdnP (2 µg)
Lane 4: CdnP (1 µg)

Representative gel image shown; actual purity may vary between batches.

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, 09/28/2022

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

PRODUCT INFORMATION



Description

Cyclic dinucleotide phosphodiesterase (CdnP) is a soluble phosphodiesterase involved in regulating cyclic dinucleotide signaling during intracellular infections of *M. tuberculosis*.¹ It is composed of an N-terminal Asp-His-His (DHH) domain and a C-terminal DHH-associated (DHHA1) domain that catalyze the hydrolysis of cyclic dinucleotides. CdnP hydrolyzes a variety of cyclic dinucleotides, including cyclic di-AMP (Item No. 17753), cyclic di-GMP (Item No. 17144), and 2'3'-cGAMP (Item No. 19887), which decreases their recognition by the stimulator of interferon genes (STING), a key mediator in the host innate immune response.²⁻⁴ Knockout of the gene encoding CdnP, *CnpB*, increases the production of IFN- β in isolated mouse bone marrow-derived macrophages (BMDMs) infected with *CnpB*^{-/-} *M. tuberculosis*, as well as decreases pulmonary and splenic bacterial burden and increases survival in *CnpB*^{-/-} *M. tuberculosis*-infected mice.⁵ Cayman's CdnP (*Mycobacterium tuberculosis* strain ATCC 25618H37Rv recombinant) protein contains N-terminal Trx- and His-tags followed by a thrombin cleavage site and an S-tag followed by an enterokinase cleavage site.

References

1. He, Q., Wang, F., Liu, S., *et al.* Structural and biochemical insight into the mechanism of Rv2837c from *Mycobacterium tuberculosis* as a c-di-NMP phosphodiesterase. *J. Biol. Chem.* **291**(27), 14386-14387 (2016).
2. Cheng, S.S., Chung, M.J., Lin, C.Y., *et al.* Phytochemicals from *Cunninghamia konishii* Hayata act as antifungal agents. *J. Agric. Food Chem.* **60**(1), 124-128 (2012).
3. Dey, R.J., Dey, B., Zheng, Y., *et al.* Inhibition of innate immune cytosolic surveillance by an *M. tuberculosis* phosphodiesterase. *Nat. Chem. Biol.* **13**(2), 210-217 (2017).
4. Jeong, H.U., Kwon, S.S., Kong, T.Y., *et al.* Inhibitory effects of cedrol, β -cedrene, and thujopsene on cytochrome P450 enzyme activities in human liver microsomes. *J. Toxicol. Environ. Health A* **77**(22-24), 1522-1532 (2014).
5. Chen, S.S., Zhang, Y., Lu, Q.L., *et al.* Preventive effects of cedrol against alopecia in cyclophosphamide-treated mice. *Environ. Toxicol. Pharmacol.* **46**, 270-276 (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