

PRODUCT INFORMATION



Calcineurin A α Isoform (human, recombinant)

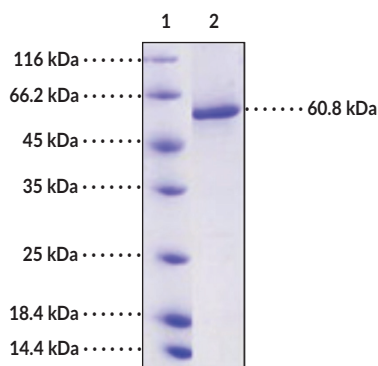
Item No. 34630

Overview and Properties

Synonyms:	Calmodulin-dependent Calcineurin A Subunit α Isoform, CALNA1, CAM-PRP Catalytic Subunit, CCN1, CNA1, PPP2B, PPP3CA, Protein Phosphatase 2B, Serine/threonine-Protein Phosphatase 2B Catalytic Subunit α Isoform
Source:	Active recombinant human N-terminal His-tagged calcineurin A expressed in insect cells
Amino Acids:	2-521 (full length)
Uniprot No.:	Q08209
Molecular Weight:	60.8 kDa
Storage:	-80°C (as supplied)
Stability:	≥ 1 year
Purity:	$\geq 94\%$ estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile 20 mM Tris, pH 8.0, with 500 mM sodium chloride and 10% glycerol
Endotoxin Testing:	< 1.0 EU/ μ g, determined by the LAL endotoxin assay
Bioactivity:	Using the Octet Red System, the affinity constant (K_d) of human PPP3CA-His bound to human PPIA-His was 0.9 nM.

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

Image



Lane 1: MW Markers
Lane 2: Calcineurin A α Isoform

SDS-PAGE Analysis of Calcineurin A α Isoform. This protein has a calculated molecular weight of 60.8 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, 10/15/2021

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

Calcineurin A is the catalytic subunit of calcineurin, a serine/threonine phosphatase with roles in T cell activation, cardiac function, learning and memory, apoptosis, and cell cycle regulation.^{1,2} Calcineurin A has three isoforms, α , β , and γ , which are encoded by *PPP3CA*, *PPP3CB*, and *PPP3CC*, respectively. Calcineurin A α is composed of an N-terminal catalytic domain, binding regions for the regulatory subunit calcineurin B (Item No. 34631) and the binding partner calmodulin, and a C-terminal autoinhibitory domain.^{3,4} It is ubiquitously expressed and associates with calcineurin B and calmodulin in a calcium-dependent manner.^{5,6} Calcineurin A α has roles in kidney development and function, as well as Tau dephosphorylation, but is dispensable for T cell development.⁶ Calcineurin A α -null mice exhibit impaired kidney function, failure to thrive, and shortened lifespan. Brain calcineurin A α levels are inversely correlated with disease severity in patients with Alzheimer's disease.⁷ Cayman's Calcineurin A α Isoform (human, recombinant) protein can be used for binding assays. This protein consists of 538 amino acids and has a calculated molecular weight of 60.8 kDa.

References

1. Rusnak, F., and Mertz, P. Calcineurin: Form and function. *Physiol. Rev.* **80(4)**, 1483-1521 (2020).
2. Mansuy, I. Calcineurin. *xPharm: The Comprehensive Pharmacology Reference*. Enna, S.J. and Bylund, D.B., editors, Elsevier (2007).
3. Gooch, J.L. An emerging role for calcineurin A α in the development and function of the kidney. *Am. J. Physiol. Renal. Physiol.* **290(4)**, F769-F776 (2006).
4. Li, H., Rao, A., and Hogan, P.G. Interaction of calcineurin with substrates and targeting proteins. *Trends Cell Biol.* **21(2)**, 91-103 (2011).
5. Nolze, A., Köhler, C., Ruhs, S., et al. Calcineurin (PPP3CB) regulates angiotensin II-dependent vascular remodelling by potentiating EGFR signalling in mice. *Acta Physiol. (Oxf.)* e13715 (2021).
6. Williams, C.R., and Gooch, J.L. Calcineurin inhibitors and immunosuppression - a tale of two isoforms. *Expert Rev. Mol. Med.* **14**, e14 (2012).
7. Karch, C.M., Jeng, A.T., and Goate, A.M. Calcium phosphatase calcineurin influences tau metabolism. *Neurobiol. Aging* **34(2)**, 374-386 (2013).

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