

Calmodulin (human), (recombinant)

Calcium-binding protein produced in *E.coli*

Citations: 14

[View Online »](#)

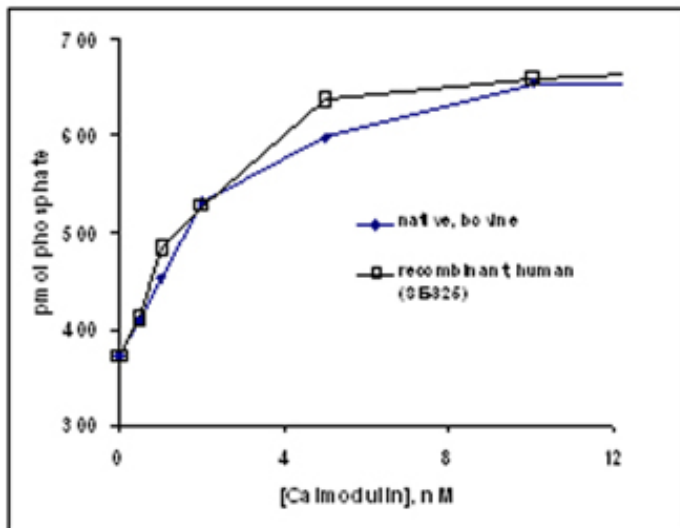
Ordering Information

[Order Online »](#)

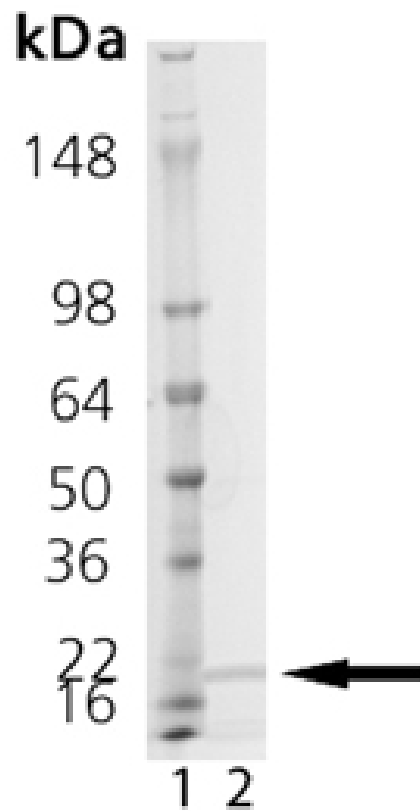
BML-SE325-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Human recombinant calmodulin (Prod. No. BML-SE325) activation of human calcineurin is essentially identical to that of purified bovine brain calmodulin (native, bovine). Initial rates of calcineurin phosphatase activity were determined as a function of calmodulin concentration with the Calcineurin Colorimetric Drug Discovery Kit (Prod. No. BML-AK804). (Note that the human and bovine calmodulin amino acid sequences are 100% identical.)



SDS-PAGE Analysis: Lane 1: MW Marker; Lane 2: 1.0 μ g of purified human recombinant calmodulin protein (Prod. No. BML-SE325).

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CaM, Phosphodiesterase 3':5'-cyclic nucleotide activator
Application Notes	Calmodulin can be used to activate calmodulin-dependent phosphodiesterase, calcineurin, CaM kinases, etc.
Biological Activity	Elicits Ca ²⁺ -dependent activation of calcineurin (Ca ²⁺ /calmodulin-dependent Ser/Thr protein phosphatase 2B, PP-2B) with the same potency and to the same maximum extent as purified native calmodulin.
CAS	73298-54-1
Formulation	Lyophilized. Salt-free.
MI	14: 1719
MW	16.8 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Reconstitution	Reconstitute in water or an appropriate buffer.
Sequence	Sequence identical to Genbank accession M27319.1.
Source	Produced in <i>E. coli</i> .
UniProt ID	P62158



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)