

Calcineurin (human), (recombinant)

Citations: 8

[View Online »](#)

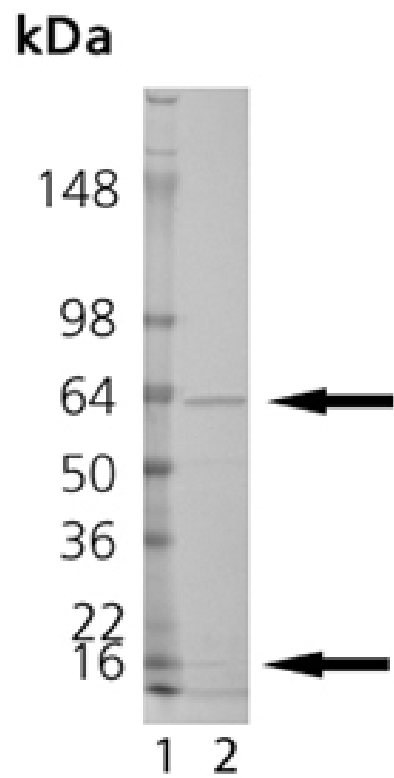
Ordering Information

[Order Online »](#)

BML-SE163-5000	5000U
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker; Lane 2: 1.0 μ g of purified human recombinant calcineurin protein (SE-163).

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Protein phosphatase 2B, PP2B
Application Notes	Useful for the study of enzyme kinetics, regulation, to dephosphorylate target substrates, and for screening inhibitors. Enzyme activity may be determined using the BIOMOL GREEN Calcineurin Assay Kit (Prod. No. BML-AK804).
Concentration	100 U/μl
Formulation	Liquid. In 1X KI-128 Assay Buffer with increased DTT concentration (50mM TRIS, pH 7.5, containing 100mM sodium chloride, 6mM MgCl ₂ , 5mM DTT, 0.025% NP-40 and 0.5mM CaCl ₂).
MW	60 kDa (calcineurin Aα) + 15 kDa (calcineurin B)
Purity	≥90% (SDS-PAGE)
Source	Produced in <i>E. coli</i> . Human calcineurin Aα (MW 60kDa) and calcineurin B (MW 15kDa) heterodimer. Both subunits are coexpressed in a construct with yeast myristoyl-CoA:protein N-myristoyltransferase. The resulting highly active calcineurin (protein phosphatase 2B) is N-myristoylated on the CaNB subunit, similar to the native protein.
Specific Activity	One U= 1pmol/min using 0.15mM RII phosphopeptide substrate (Prod. No. BML-P160) in the presence of calcium/calmodulin at 30°C. In the absence of calmodulin, the activity is <1%.
UniProt ID	Q08209 (subunit Aα), P63098 (subunit Bα)



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)