

Calcineurin cellular activity assay kit

Provides everything needed to assay calcineurin cellular activity, including recombinant calcineurin control

A complete colorimetric assay kit for measuring cellular calcineurin (PP2B) phosphatase activity. It employs a convenient 96-well plate format with all reagents necessary for measuring calcineurin (PP2B) phosphatase activity in tissue/cellular extracts PLUS human recombinant calcineurin is included as a positive control. The RII phosphopeptide substrate, supplied with this kit is the most efficient and selective peptide known for calcineurin. The detection of free-phosphate released is based on the classic malachite green assay.

Citations: 68

[View Online »](#)

Ordering Information

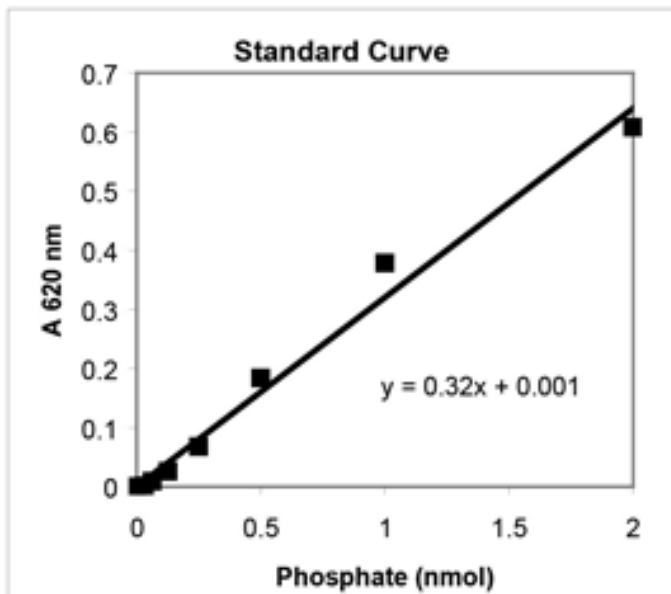
[Order Online »](#)

BML-AK816-0001	96 wells
-----------------------	----------

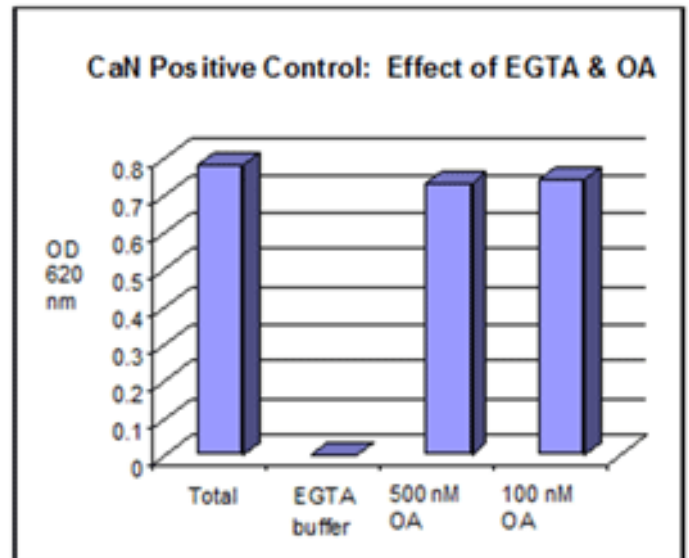
Manuals, SDS & CofA

[View Online »](#)

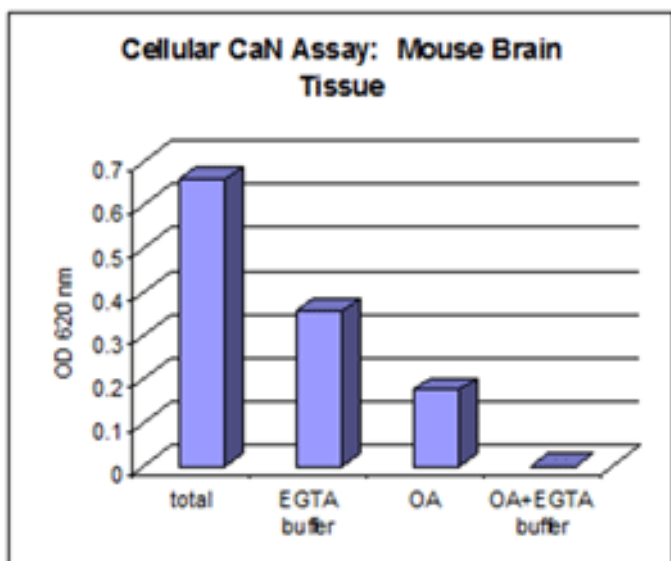
- Safe and non-radioactive assay
- Convenient 1-step detection
- Excellent sensitivity



BIOMOL GREEN™ Phosphate Standard Curve



CaN Positive Control. Purified calcineurin (BML-SE163-9090) positive control. The CaN was incubated 1 hr at 30°C under various buffer conditions. The results demonstrate that CaN activity is inhibited by the EGTA buffer, but not inhibited by 100nM or 500nM concentrations of okadaic acid



Cellular CaN Assay Using Mouse Brain Phosphatase activity from a freshly prepared mouse brain extract. Prior to gel filtration the extract was diluted 1:1 in lysis buffer containing protease inhibitors. After gel filtration and prior to the phosphatase assay, the extract was diluted 1/25 in lysis buffer. Well contents were as in manual. The reaction was incubated for 30 min at 30°C.

Handling & Storage

Use/Stability	Store all components except the microtiter plate at 80°C for the highest stability. The calcineurin enzyme component SE163-9090 must be handled particularly carefully in order to retain maximal enzymatic activity. Thaw it quickly in a RT water bath or by rubbing between fingers, then immediately store on an ice bath. The remaining unused enzyme should be quickly refrozen by placing at -80°C. To minimize the number of freeze/thaw cycles, aliquot the calcineurin into separate tubes and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	Colorimetric detection
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	Calcineurin enzyme, Calmodulin (human, recombinant), Substrate (R11 phosphopeptide), 2x assay buffer, EGTA Buffer, Lysis Buffer, Protease inhibitor cocktail, Biomol Green™ Reagent, Phosphate standard, Okadaic acid, Desalting column and resin, 1/2-Volume microplate



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)