

Calcineurin Phosphatase assay kit

Designed for inhibitor screening or kinetic assay of recombinant Calcineurin

This colorimetric kit is designed for inhibitor screening or kinetic assay of recombinant Calcineurin. Activity is determined using the RII phosphopeptide substrate, the most efficient and selective peptide known for calcineurin, and detection of free phosphate released is based on the classic malachite green assay.

Citations: 17

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AK804-0001	96 wells
----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Safe and non-radioactive assay
- Convenient 1-step detection, no mixing needed
- High throughput with 96 well plate format

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Activity assay, Colorimetric detection, HTS

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Recombinant calcineurin, RII phosphopeptide substrate, calmodulin, assay buffer, BIOMOL[®] GREEN detection reagent, phosphate standard, microplate and detailed instructions.



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)