

[pSer95, Ala97]-RII (81-99)

Calcineurin substrate

This sequence, DLDVPIPGRFDRRVpSVAAE is derived from cAMP-dependent protein kinase regulatory subunit type II [PKA; [Ala-97]-RII (81-99)] with a phosphoserine at Ser⁹⁵. This peptide can be used directly as a calcineurin (protein phosphatase 2B, PP2B) substrate. Phosphate release can be quantified using the (BML-AK111) detection reagent.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P160-0500	0.5mg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solution in water is highly stable at -20°C
Handling	Protect from moisture
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	RII Phosphopeptide, [pSer95, Ala97]-cAMP-dependent protein kinase regulatory subunit type II (81-99)
Appearance	Solid.
Formulation	Lyophilized. Trifluoroacetate salt.
Identity	Determined by MS.
MW	2192.0
Purity	≥95% (HPLC)
Sequence	Asp-Leu-Asp-Val-Pro-Ile-Pro-Gly-Arg-Phe-Asp-Arg-Arg-Val- pSer -Val-Ala-Ala-Glu
Solubility	Soluble in water.
Source	Synthetic.

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)