

CaM kinase IV substrate

This sequence, KSDGGVKKRKSSSS, is from Ca^{2+} /calmodulin-dependent protein kinase II- γ (345-358). Substrate is specific for CaMKIV ($K_m = 8 \mu\text{M}$) and is a poor substrate for type II. Useful with purified enzyme and crude extracts using phosphocellulose assay.

Ordering Information

[Order Online »](#)

BML-P100-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Ca²⁺/calmodulin-dependent protein kinase II γ (345-358), Peptide γ , Ca²⁺/calmodulin-dependent protein kinase IV substrate, CAMKIV substrate

MW 1450.8

Sequence Lys-Ser-Asp-Gly-Gly-Val-Lys-Lys-Arg-Lys-Ser-Ser-Ser-Ser

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-
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)