

[Ala⁹⁷]-RII (81-99)

PKA substrate

This sequence, DLDVPIPIGRFDRRVSVAEE, is from cAMP-dependent protein kinase regulatory subunit type II (PKA). The phosphorylated peptide (BML-P160) can be used as a substrate for calcineurin ($K_m = 25 \mu\text{M}$, $V_{\text{max}} = 6 \mu\text{mol}/\text{min}/\text{mg}$).

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	[Ala97]-cAMP-dependent protein kinase regulatory subunit type II (81-99)
MW	2113.0
Purity	≥97%
Sequence	Asp-Leu-Asp-Val-Pro-Ile-Pro-Gly-Arg-Phe-Asp-Arg-Arg-Val-Ser-Val-Ala-Ala-Glu

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com