

BPDEtide

Substrate for PKG and PKA

Sequence derived from bovine lung cGMP specific phosphodiesterase (cG-BPDE). Specific substrate for PKG, $K_m=68 \mu\text{M}$; K_m for PKA=320 μM , with no phosphorylation detected for CaMKII or PKC.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P112-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	cGMP-dependent protein kinase (PKG) substrate
Appearance	White to off-white solid.
Formula	$C_{68}H_{115}N_{23}O_{20}$
Formulation	Lyophilized.
MW	1574.8
Purity	≥97% (HPLC)
Sequence	H-Arg-Lys-Ile-Ser-Ala-Ser-Glu-Phe-Asp-Arg-Pro-Leu-Arg-OH
Solubility	Soluble in water.
Source	Synthetic.

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-fr@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