

MARCKS protein (bovine brain) (151- 175)

Excellent PKC substrate

Excellent substrate for protein kinase C (PKC) ($K_m = 20\text{nM}$ for purified rat brain PKC). The peptide consists of residues 151-175 of the myristoylated alanine-rich C-kinase substrate (MARCKS) protein phosphorylation site. Protein kinase C isotypes (including PKC ϵ) phosphorylate serines 8, 12 and 19 of this peptide. It is a very weak substrate for cAMP- and cGMP-dependent protein kinase and CaM kinases I, II and III.

Citations: 2

[View Online »](#)

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	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white powder.
CAS	125267-21-2
Formula	$C_{147}H_{243}N_{41}O_{31}$
Formulation	Lyophilized
MW	3081.8
Purity	≥98%
Sequence	H-Lys-Lys-Lys-Lys-Lys-Arg-Phe-Ser-Phe-Lys-Lys-Ser-Phe-Lys-Leu-Ser-Gly-Phe-Ser-Phe-Lys-Lys-Asn-Lys-Lys-OH
Solubility	Soluble in water.

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