

# 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 $\epsilon$ ) 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 »](#)

## Ordering Information

[Order Online »](#)

ALX-162-004-M001	1mg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Appearance</b>	White to off-white powder.
<b>CAS</b>	125267-21-2
<b>Formula</b>	$C_{147}H_{243}N_{41}O_{31}$
<b>Formulation</b>	Lyophilized
<b>MW</b>	3081.8
<b>Purity</b>	≥98%
<b>Sequence</b>	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
<b>Solubility</b>	Soluble in water.



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)