

# Kemptide

## PKA substrate

Phosphate acceptor peptide. Substrate for Ser/Thr protein kinases.  $K_m = 4.7 \mu\text{M}$  for cAMP-dependent protein kinase (PKA). The sequence defines the phosphorylation site of porcine liver pyruvate kinase.

Citations: 8

[View Online »](#)

## Ordering Information

[Order Online »](#)

|               |     |
|---------------|-----|
| BML-P107-0001 | 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

**CAS** 65189-71-1

**Formula**  $C_{32}H_{61}N_{13}O_9$

**Formulation** Lyophilized.

**MW** 771.9

**Purity** ≥97% (HPLC)

**Sequence** H-Leu-Arg-Arg-Ala-Ser-Leu-Gly-OH

**Solubility** Soluble in water.

**Technical Info / Product Notes** Replacement for **ADI-SPK-101**

**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)