

# MLCK substrate (skeletal and smooth muscle)

This sequence, KKRAARATSDVFA, is chicken gizzard smooth muscle myosin P-light chain (MLC 11-23) with Pro to Ala<sup>14</sup> and Gln to Ala<sup>15</sup> substitutions. Contains specificity determinants for MLC phosphorylation (target is Ser<sup>19</sup>) and is an excellent myosin light chain kinase (MLCK) substrate.  $K_m = 8.6 \mu\text{M}$  for rabbit skeletal muscle MLCK,  $7.5 \mu\text{M}$  for chicken gizzard smooth muscle MLCK.

## Ordering Information

[Order Online »](#)

BML-P114-0001	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	Myosin light chain kinase substrate, Lys-Lys-Arg-Ala-Ala-Arg-Ala-Thr-Ser-Asn-Val-Phe-Ala-NH <sub>2</sub>
Appearance	White to off-white solid.
Formulation	Lyophilized.
Identity	Determined by MS.
MW	1418.0
Purity	≥97% (HPLC)
Sequence	Lys-Lys-Arg-Ala-Ala-Arg-Ala-Thr-Ser-Asn-Val-Phe-Ala-NH <sub>2</sub>
Solubility	Soluble in water.
Source	Synthetic.

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-  
fr@enzolifesciences.com](mailto:info-fr@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)