

MLCK substrate (skeletal and smooth muscle)

This sequence, KKRAARATSDVFA, is chicken gizzard smooth muscle myosin P-light chain (MLC 11-23) with Pro to Ala¹⁴ and Gln to Ala¹⁵ substitutions. Contains specificity determinants for MLC phosphorylation (target is Ser¹⁹) 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 ₂
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 ₂
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-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