

MLCK (342-352) amide

MLCK inhibitor

This sequence, KRRWKKNFIAV, is from skeletal muscle myosin light chain kinase (MLCK) (342-352) pseudosubstrate site. Inhibits MLCK ($K_i = 1 \mu\text{M}$ with MLC peptide substrate). Possesses only slight calmodulin binding activity.

Ordering Information

[Order Online »](#)

BML-P209-0001	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

Alternative Name Myosin light chain kinase (342-352) amide, MLCK inhibitor

MW 1444.9

Purity ≥97%

Sequence Lys-Arg-Arg-Trp-Lys-Lys-Asn-Phe-Ile-Ala-Val-NH₂

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)