

Casein kinase II β

(198-215)

CDK1 substrate

This sequence, QLQLQAASDFKSPVKTI, contains a CKII autophosphorylation site plus a CDK1 (p34^{Cdc2}) consensus phosphorylation site Ser²⁰⁹. Used as a CDK1 substrate (0.25 mM) and also a competitive inhibitor with IC₅₀=50 μ M for inhibition of apoptosis induced by perforin and fragmentin-2 in YAC-1 lymphoma cells (peptide A).

Ordering Information

[Order Online »](#)

BML-P103-0001	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

Appearance	Lyophilized solid.
MW	2028.1
Sequence	Gln-Leu-Gln-Leu-Gln-Ala-Ala-Ser-Asn-Phe-Lys-Ser-Pro-Val-Lys-Thr-Ile-Arg

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