

Casein kinase II (human), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE124-0010	10KU
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CK II
Formulation	Liquid. In 20mM TRIS-HCl, pH 7.5, containing 0.35M sodium chloride, 1 mM EDTA, 2 mM DTT and 0.1% Triton X-100.
MW	44 and 26 kDa subunits
Purity Detail	Free from proteases and phosphatases.
Source	Produced in <i>E. coli</i> . Consists of 44kDa α catalytic and 26kDa β regulatory subunits as an $\alpha_2\beta_2$ tetramer.
Specific Activity	One unit will transfer 1 pmol phosphate from ATP to peptide substrate per minute at 30°C.

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