

PKA α (catalytic domain) (mouse), (recombinant)

Ordering Information

[Order Online »](#)

BML-SE192-5000	5000U
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	cAMP-dependent protein kinase α , Protein kinase A α
Application Notes	Useful for phosphorylation of target regulatory enzymes and other substrates of PKA.
Concentration	~0.5mg/ml
Formulation	Liquid. In 50% glycerol buffer.
MW	38 kDa
Purity Detail	Contains no detectable protease or phosphatase activity.
Source	Produced in <i>E. coli</i> .
Specific Activity	One unit will add 1 nmol phosphate to kemptide peptide substrate (P-107) per hour at pH 7.5, 30°C. Regulatory subunit is not present and therefore activity is not dependent on cAMP.
UniProt ID	P05132

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)