

MKP-3 (human), (recombinant)

MKP-3 is a member of a group of dual specificity phosphatases which negatively regulate MAP kinases, dephosphorylating both the phosphotyrosine and phosphothreonine in the pTXpY activation loop motif. Transcriptional suppression of MKP-3 expression has been implicated in the development of pancreatic cancer. MKP-3 is located in the nucleus where it may act to anchor Erk 1/2.

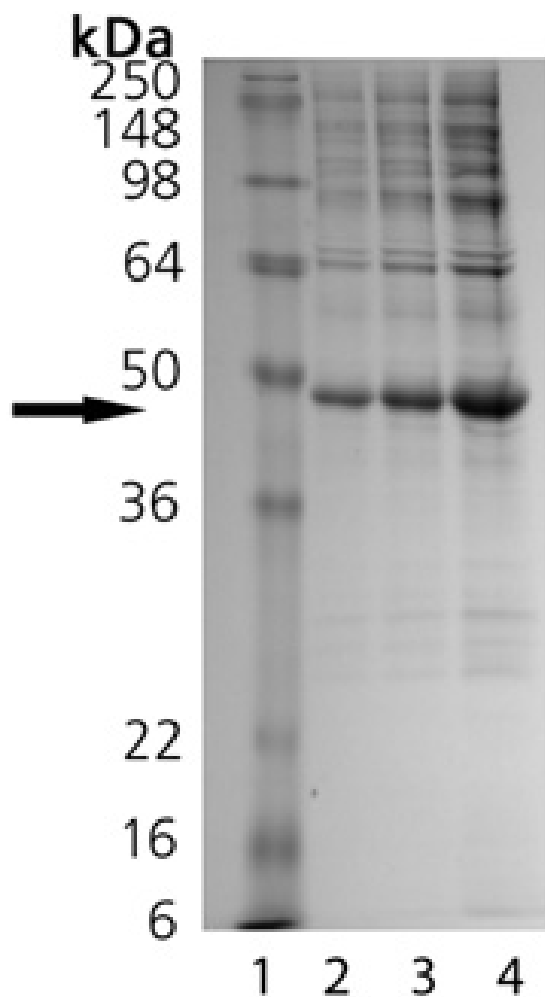
Ordering Information

[Order Online »](#)

BML-SE454-0100	100µg
-----------------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis. Lane 1: MW Marker, Lane 2: 1 µg, Lane 3: 2 µg, Lane 4: 5 µg MKP-3.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name

MAP kinase phosphatase 3, DUSP6, Dual specificity protein phosphatase 6, PYST1

Formulation

Liquid. In 20mM TRIS/HCl, pH 8.0, 200mM NaCl, 5mM DTT, 0.1 % Tween-20, 10% glycerol.

Gene/Protein Identifier

NM_001946 (RefSeq)

MW

42 kDa

Purity Detail

Partially purified by single-step affinity chromatography and gel filtration.

Source

Produced in Sf9 insect cells. Recombinant full length MKP-3. Produced in a baculovirus expression system.

Specific Activity

≥10 U/mg assayed by 3-O-methyl fluorescein phosphate (OMFP) hydrolysis at pH 8.2, 30°C. One unit is equal to 1 nmole phosphate hydrolyzed from OMFP per minute.

UniProt ID

Q16828

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-
fr@enzolifesciences.com](mailto:info-fr@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)