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.

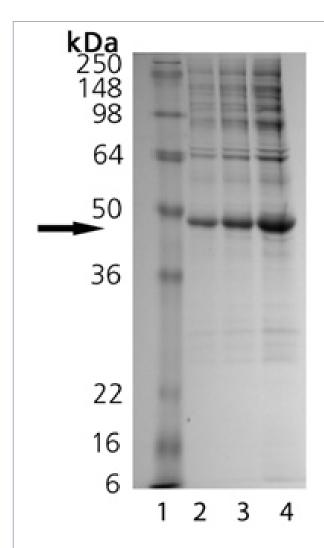
Ordering Information

Order Online »

BML-SE454-0100 100μg

Manuals, SDS & CofA

View Online »



SDS-PAGE Analysis. Lane 1: MW Marker, Lane 2: 1 μ g, Lane 3: 2 μ g, Lane 4: 5 μ g 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

