

# MKP-5 (human), (recombinant) (His- tag)

MKP-5 (DUSP10) 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. MKP5 has a principal function in both innate and adaptive immune responses.

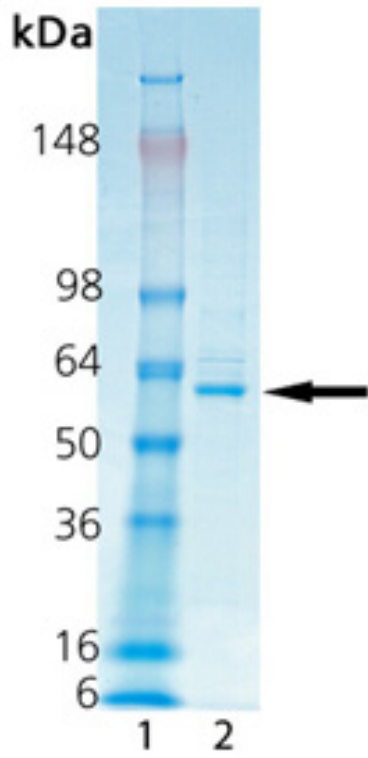
## Ordering Information

[Order Online »](#)

BML-SE467-0100	100µg
----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis. Lane 1: MW Marker, Lane 2: 1.0 $\mu$ g of purified MKP-5 (human) (recombinant) (His-tag).

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	MAP kinase phosphatase 5, DUSP10, Dual specificity protein phosphatase 10
<b>Formulation</b>	Liquid. In 20mM TRIS/HCl, pH 8.0, 200mM NaCl, 5mM DTT, 0.1 % Tween-20, 10% glycerol.
<b>Gene/Protein Identifier</b>	NM_007207 (RefSeq)
<b>MW</b>	53 kDa
<b>Purity Detail</b>	Partially purified by single-step affinity chromatography and gel filtration.
<b>Source</b>	Produced in Sf9 insect cells. Recombinant human MKP-5 fused at the N-terminus to a His-tag. Produced in a baculovirus expression system.
<b>Specific Activity</b>	≥50 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.
<b>UniProt ID</b>	Q9Y6W6

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)