

VHR dual specificity phosphatase (human), (recombinant)

Dual specificity phosphatases (DUSPs) are members of the superfamily of protein tyrosine phosphatases. These phosphatases inactivate their target kinases by dephosphorylating both the phosphoserine/threonine and phosphotyrosine residues. They negatively regulate members of the mitogen-activated protein (MAP) kinase superfamily (MAPK/ERK, SAPK/JNK, p38), which are associated with cellular proliferation and differentiation.

Citations: 2

[View Online »](#)

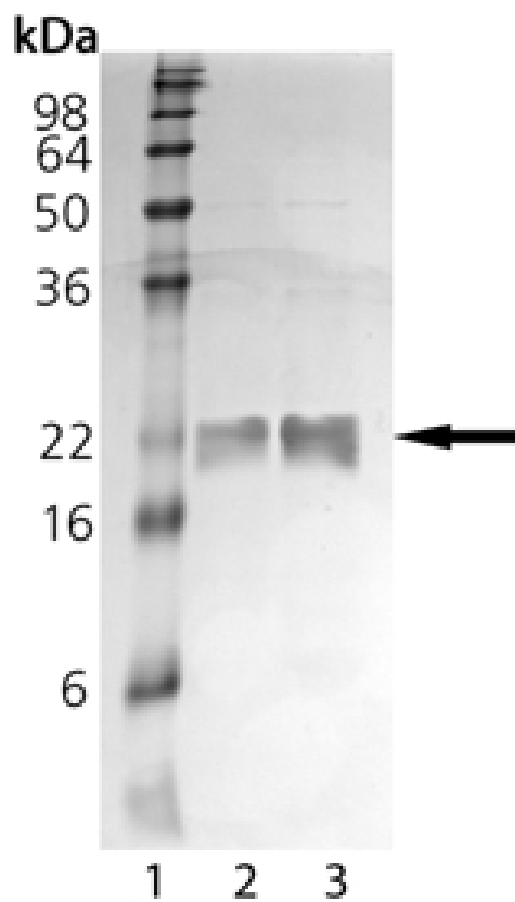
Ordering Information

[Order Online »](#)

BML-SE333-0050	50µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 μ g, Lane 3: 2 μ g VHR dual-specificity phosphatase

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 Dual specificity protein phosphatase 3, DUSP3

Application Notes Useful for the study of VHR dual phosphatase kinetics and substrate specificity, and to dephosphorylate pSer/Thr and pTyr phosphorylated protein substrates.

Formulation Liquid. In 50mM citrate, pH 6.0, containing 0.1M sodium chloride, 1mM EDTA and 1mM DTT.

MW 20.4 kDa

Purity ≥95% (SDS-PAGE)

Source Produced in *E. coli*.

Specific Activity 11U/mg. One unit will hydrolyze 1μmol p-nitrophenyl phosphate (pNPP) per minute at pH 6.0 and 30°C. Reaction conditions are 50mM citrate, pH 6.0, 0.1M NaCl, 1mM EDTA and 1mM DTT, 10mM pNPP substrate. An increase in signal of 3 to 4 fold can be achieved by addition of base to a pH of approximately 9.0 at the end of an assay with pNPP.

UniProt ID P51452

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