

VHR dual specificity phosphatase (human), (recombinant)

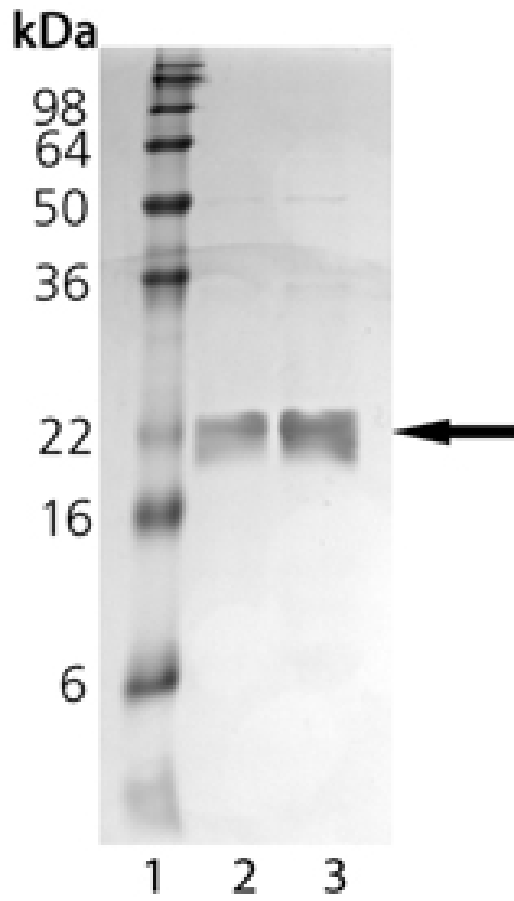
Dual specificity phosphatases (DUSPs) are members of the superfamily of protein tyrosine phosphatases. These phosphatases inactivate their target kinases by dephosphorylating both the phosphoserine/threonine and phosphotyrosine residues. They negatively regulate members of the mitogen-activated protein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

Citations: 2

[View Online »](#)

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 <i>E. coli</i> .
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](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)