

# PRL-1 (human), (recombinant) (His- tag)

PRL phosphatases comprise a class of small oncogenic phosphatases that are prenylated at their carboxyl-termini. PRL-1 is overexpressed during cell proliferation in the liver and during differentiation of epithelial cells in the digestive tract. It augments MMP-2 and MMP-9 expression, which in turn induce cell migration and invasion.

## Ordering Information

[Order Online »](#)

BML-SE556-0050	50µg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Dilution of the enzyme followed by refreezing may lead to loss of activity.
Long Term Storage	-80°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Phosphatase of regenerating liver 1, Protein tyrosine phosphatase type IVA 1
Application Notes	Useful for studies of enzyme kinetics and regulation, dephosphorylation of target substrates, and inhibitor screening.
Formulation	Liquid. In 40mM TRIS-HCl, pH 8.0, containing 110mM sodium chloride, potassium chloride, 200mM imidazole, 20% glycerol and 3mM DTT.
MW	~19.8kDa
Purity	≥75% (SDS-PAGE)
Source	Produced in <i>E. coli</i> . The catalytic domain of PRL-1 (aa 2-173) is fused at the N-terminus to a His-tag.
Specific Activity	≥0.168U/min/μg. One unit will hydrolyze 1pmol <i>para</i> -nitrophenyl phosphate (PNPP) per minute at 37°C. Assay buffer: 50mM TRIS, pH 7.4, containing 150mM sodium chloride, 5mM dithiothreitol and 12.5mM PNPP.
UniProt ID	Q93096

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)