

# LAR (human), (recombinant)

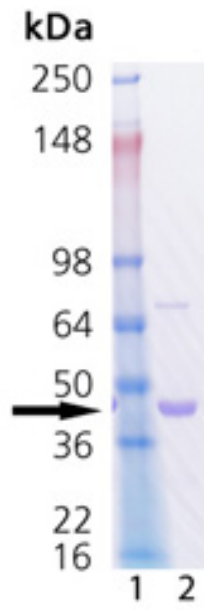
Removes phosphate specifically from tyrosine residues in proteins. Rate of dephosphorylation depends on LAR substrate preference.

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 $\mu$ g LAR protein tyrosine phosphatase (human), (recombinant)

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	Leukocyte common antigen-related protein tyrosine phosphatase
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.0, containing 100mM sodium chloride, 2mM EDTA, 5mM dithiothreitol, 0.01% Brij 35, and 30% glycerol.
MW	37 kDa
Purity	≥90% (SDS-PAGE)
Purity Detail	Contains no detectable protease, DNase, RNase, or Ser/Thr phosphatase activity.
Source	Produced in <i>E. coli</i> . Soluble catalytic LAR-D1 domain (350 aa).
Specific Activity	≥3 U/μg. One unit will hydrolyze 1 nmol p-nitrophenyl phosphate per minute at pH 7.0 and 30°C. 20 U can dephosphorylate >95% of phosphotyrosine in 1 nmol protein in 30 minutes.
UniProt ID	P10586

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)