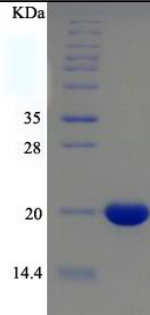


Recombinant Human ACP1 Protein Data Sheet

Catalog #	hRP-C0027-EF011
Size	25 µg
Protein Name	Human low molecular weight phosphotyrosine protein phosphatase
Protein Symbol	ACP1
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	NM_004300.1
Uniprot Accession #	P24666
Description	Acid phosphatase1 belongs to the phosphotyrosine protein phosphatase family of proteins. It functions as an acid phosphatase and a protein tyrosine phosphatase by hydrolyzing protein tyrosine phosphate to protein tyrosine and orthophosphate. This enzyme also hydrolyzes orthophosphoric monoesters to alcohol and orthophosphate.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	173aa(including fusion tag)
Molecular Weight	20.0kDa(including fusion tag)
pI	7.2
Activity	NA
Storage	Storage buffer: 20mM Tris.Cl, 50mM NaCl, 50% Glycerol, pH9.0. Store at -80°C and avoid repeated freeze-thaw cycles.
	
Reference:	• Low-molecular-weight protein tyrosine phosphatase and human disease: in search of biochemical mechanisms • From immune response to cancer: a spot on the low molecular weight protein tyrosine phosphatase • The low-molecular-weight phosphotyrosine phosphatase is a negative regulator of FcγRIIA-mediated cell activation • Regulation of the EphA2 kinase by the low molecular weight tyrosine phosphatase induces transformation • Low molecular weight protein tyrosine phosphatase (LMW-PTP) as a diagnostic and therapeutic target

GeneCopoeia™
Expressway to Discovery

GeneCopoeia Inc.

9620 Medical Center Drive, Suite 101

Rockville, MD 20850

USA

Phone: 301-762-0888

Toll free: 1-301-762-0888

Fax: 301-762-3888

Web: www.genecopoeia.com

Inquiry: inquiry@genecopoeia.com

Technical Support: support@genecopoeia.com