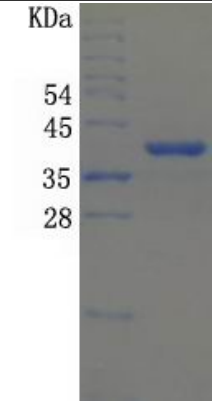


Recombinant Human ARG1 Protein Data Sheet

Catalog #	hRP-C0062-EF012
Size	100 µg
Protein Name	Human arginase-1
Protein Symbol	ARG1
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	NM_000045.2
Uniprot Accession #	P05089
Description	Arginase catalyzes the hydrolysis of arginine to ornithine and urea. At least two isoforms of mammalian arginase exist(types I and II) which differ in their tissue distribution, subcellular localization, immunologic cross reactivity and physiologic function. The type I isoform encoded by this gene, is a cytosolic enzyme and expressed predominantly in the liver as a component of the urea cycle. Inherited deficiency of this enzyme results in argininemia, an autosomal recessive disorder characterized by hyperammonemia.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	337aa(including fusion tag)
Molecular Weight	36.6kDa(including fusion tag)
pI	7.5
Activity	NA
Storage	Storage buffer: 1XPBS. Store at -80°C and avoid repeated freeze-thaw cycles.
	 <p>The image shows an SDS-PAGE gel with a single prominent band at approximately 36.6 kDa. The molecular weight markers on the left are 54, 45, 35, and 28 kDa. The band is located between the 35 and 45 kDa markers, consistent with the stated molecular weight of 36.6 kDa including the fusion tag.</p>

Reference:	<ul style="list-style-type: none">• Vascular dysfunction in retinopathy-an emerging role for arginase• Structure and function of arginases• Arginase I deficiency: severe infantile presentation with hyperammonemia: more common than reported• Arginase-1: a new immunohistochemical marker of hepatocytes and hepatocellular neoplasms• Functionally important role for arginase 1 in the airway hyperresponsiveness of asthma
------------	---



GeneCopoeia Inc.

9620 Medical Center Drive, Suite 101
Rockville, MD 20850
USA

Phone: 301-762-0888

Toll free: 1-866-360-9531

Fax: 301-762-3888

Web: www.genecopoeia.com

Inquiry: inquiry@genecopoeia.com

Technical Support: support@genecopoeia.com