Recombinant Human ARG1 Protein Data Sheet

Catalog #	hRP-C0062-EF012				
Size	100 μg				
Protein Name	Human arginase-1				
Protein Symbol	ARG1				
Original Source	Homo sapiens				
Expression System	E.coli				
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 inthe 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.				
	KDa 54 45 35 28				
	Vascular dysfunction in retinopathy-an emerging role for arginase				
Reference:	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: 1-301-762-0888 Fax: 301-762-3888

Web: www.genecopoeia.com
Inquiry: inquiry@genecopoeia.com

 $Technical \ Support: \underline{support@genecopoeia.com}$