Recombinant Human AKR1A1 Protein Data Sheet

Catalog #	hRP-C0052-EF012
Size	100 µg
Protein Name	Human aldo-keto reductase family 1 member A1
Protein Symbol	AKR1A1
Original Source	Homo sapiens
Expression System	E.coli
GenBank Accession #	NM 006066.1
Uniprot Accession #	P14550
Description	This gene encodes a member of the aldo/keto reductase superfamily, which consists of more than 40 known enzymes and proteins. This member, also known as aldehyde reductase, is involved in the reduction of biogenic and xenobiotic aldehydes and is present in virtually every tissue. Multiple alternatively spliced transcript variants of this gene exist, all encoding the same protein.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	340aa(including fusion tag)
Molecular Weight	38.5 kDa (including fusion tag)
pI	7.1
Activity	NA
Storage	Storage buffer: 20mM Tris.Cl, 100mM NaCl, 0.4M Urea, 50% Glycerol, pH8.0. Store at -80°C and avoid repeated freeze-thaw cycles.
	67 54 45 35 28
Reference:	 Mechanism of human aldehyde reductase: characterization of the active site pocket Quantitative determination of human aldose reductase by enzyme-linked immunosorbent assay: Immunoassay of human aldose reductase Aldose Reductase in Glucose Toxicity: A Potential Target for the Prevention of Diabetic Complications Enzyme immunoassay for erythrocyte aldose reductase



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