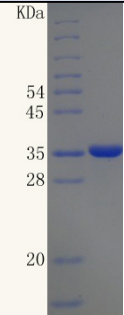


Recombinant Human AKR1B1 Protein Data Sheet

Catalog #	hRP-C0237-EF011
Size	25 µg
Protein Name	Human aldo-keto reductase family 1 member B1
Protein Symbol	AKR1B1
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	NM_001628.1
Uniprot Accession #	P15121
Description	This gene encodes a member of the aldo/keto reductase superfamily, which consists of more than 40 known enzymes and proteins. This member catalyzes the reduction of a number of aldehydes, including the aldehyde form of glucose, and is thereby implicated in the development of diabetic complications by catalyzing the reduction of glucose to sorbitol. Multiple pseudogenes have been identified for this gene. The nomenclature system used by the HUGO Gene Nomenclature Committee to define human aldo-keto reductase family members is known to differ from that used by the Mouse Genome Informatics database.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	331aa(including fusion tag)
Molecular Weight	37.8kDa(including fusion tag)
pI	7.2
Activity	NA
Storage	Storage buffer: 20mM Tris.Cl, 50mM NaCl, 50% Glycerol, pH8.0. Store at -80°C and avoid repeated freeze-thaw cycles.
	
Reference:	• Targeting aldose reductase for the treatment of cancer. • Understanding the role of aldose reductase in ocular inflammation • Aldose reductase and cardiovascular diseases, creating human-like diabetic complications in an experimental model. • The crystal structure of the aldose reductase.NADPH binary complex. • Aldose reductase: a novel therapeutic target for inflammatory pathologies. • Aldose reductase in diabetic microvascular complications • Quantitative determination of human aldosereductase by enzyme-linked immunosorbent assay: Immunoassay of human aldosereductase

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