## Recombinant Human AKR1A1 Protein Data Sheet

Catalog #	hRP-C0052-EF011				
Size	25 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:	<ul> <li>Mechanism of human aldehyde reductase: characterization of the active site pocket</li> <li>Quantitative determination of human aldose reductase by enzyme-linked immunosorbent assay: Immunoassay of human aldose reductase</li> <li>Aldose Reductase in Glucose Toxicity: A Potential Target for the Prevention of Diabetic Complications</li> <li>Enzyme immunoassay for erythrocyte aldose reductase</li> </ul>				



## 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: <a href="mailto:www.genecopoeia.com">www.genecopoeia.com</a>
Inquiry: <a href="mailto:inquiry@genecopoeia.com">inquiry@genecopoeia.com</a>

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