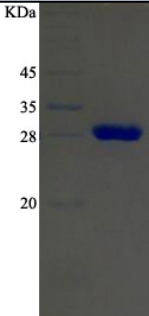


Recombinant Human AK2 Protein Data Sheet

Catalog #	hRP-X0417-EF011
Size	25 µg
Protein Name	Human adenylate kinase 2, mitochondrial
Protein Symbol	AK2
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	NM_001625.1
Uniprot Accession #	P54819
Description	Three isozymes of adenylate kinase, namely 1, 2, and 3, have been identified in vertebrates. They are encoded by transcript variants derived from the same gene. Expression of these isozymes is tissue-specific and developmentally regulated. Isozyme 2 is localized in the mitochondrial intermembrane space and may play a role in apoptosis.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	254aa(including fusion tag)
Molecular Weight	28.4kDa(including fusion tag)
pI	8.0
Activity	NA
Storage	Storage buffer: 20mM Tris.Cl, 50mM NaCl, 20% Glycerol, pH7.0. Store at -80°C and avoid repeated freeze-thaw cycles.
	
Reference:	• Adenylate kinase and AMP signaling networks: Metabolic monitoring, signal communication and body energy sensing • Human adenylate kinase 2 deficiency causes a profound hematopoietic defect associated with sensorineural deafness • Adenylate kinase 2, a mitochondrial enzyme • AK2 activates a novel apoptotic pathway through formation of a complex with FADD and caspase-10 • Dynamics of nucleotide metabolism as a supporter of life phenomena • Cloning and Expression of Human Adenylate Kinase 2 Isozymes :Differential Expression of Adenylate Kinase 1 and 2 in Human Muscle Tissues

GeneCopoeia™
Expressway to Discovery

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