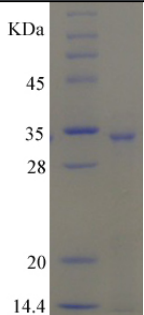


Recombinant Human IL1A Protein Data Sheet

Catalog #	hRP-A0179-EF011
Size	25 µg
Protein Name	Human mRNA for interleukin-1 precursor
Protein Symbol	IL1A
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	X02851.1
Uniprot Accession #	P01583
Description	Produced by activated macrophages, IL-1 stimulates thymocyte proliferation by inducing IL-2 release, B-cell maturation and proliferation, and fibroblast growth factor activity. IL-1 proteins are involved in the inflammatory response, being identified as endogenous pyrogens, and are reported to stimulate the release of prostaglandin and collagenase from synovial cells.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	286aa
Molecular Weight	32.5kDa
pI	5.7
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.
	
Reference:	• Structure of interleukin 1 alpha at 2.7-A resolution. • Association of IL1A gene polymorphisms with susceptibility to and severity of systemic sclerosis in the Japanese population

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