## 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	Homo sapiens
Expression System	E.coli
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.
	KDa 45 35 28 20 14.4
Reference:	<ul> <li><u>Structure of interleukin 1 alpha at 2.7-A resolution.</u></li> <li><u>Association of IL1A gene polymorphisms with susceptibility to and severity of systemic sclerosis in the Japanese population</u></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: www.genecopoeia.com Inquiry: inquiry@genecopoeia.com Technical Support: support@genecopoeia.com