Recombinant Human IL1B Protein Data Sheet

Catalog #	hRP-A0180-ES131
Size	25 μg
Protein Name	Human interleukin-1 beta
Protein Symbol	IL1B
Original Source	Homo sapiens
Expression System	E.coli
GenBank Accession #	X02532.1
Uniprot Accession #	P01584
Description	The protein encoded by this gene is a member of the interleukin 1 cytokine family. This cytokine is produced by activated macrophages as a proprotein, which is proteolytically processed to its active form by caspase 1 (CASP1/ICE). This cytokine is an important mediator of the inflammatory response, and is involved in a variety of cellular activities, including cell proliferation, differentiation, and apoptosis. The induction of cyclooxygenase-2 (PTGS2/COX2) by this cytokine in the central nervous system (CNS) is found to contribute to inflammatory pain hypersensitivity. This gene and eight other interleukin 1 family genes form a cytokine gene cluster on chromosome 2.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	Trx-His-SUMO
Peptide Length	493aa(including fusion tag)
Molecular Weight	55.6kDa(including fusion tag)
pI	4.8
Activity	NA
Storage	Storage buffer: 1XPBS. Store at -80°C and avoid repeated freeze-thaw cycles.
	kDa 60 50 40 30 25 20 • Interleukin-1 beta targeted therapy for type 2 diabetes
Reference:	• Interleukin-1 beta targeted therapy for type 2 diabetes
	• Interleukin-1 beta and interleukin-18: regulation and activity in local inflammation
	Levels of IL-1 beta control stimulatory/inhibitory growth of cancer cells
	Association study between interleukin 1 beta gene and epileptic disorders: a HuGe review and meta-analysis
	• Structural and functional comparison of cytokine interleukin-1 beta from chicken and human
	Role of interleukin 1 beta in esophageal squamous cell carcinoma



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

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