Recombinant Human IL6 Protein Data Sheet

Catalog #	hRP-A0301-EF011
Size	25 μg
Protein Name	Human interleukin 6
Protein Symbol	IL6
Original Source	Homo sapiens
Expression System	E.coli
GenBank Accession #	M14584.1
Uniprot Accession #	P05231
Description	It also known as interferon-beta 2 and B cell stimulatory factor-2 (BSF-2), is a multifunctional alpha-helical cytokine that plays important roles in host defense, acute phase reactions, immune responses, and hematopoiesis. IL-6 is expressed by a variety of normal and transformed cells including T cells, B cells, monocytes/macrophages, fibroblasts, hepatocytes, keratinocytes, astrocytes, vascular endothelial cells, and various tumor cells.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	227aa(including fusion tag)
Molecular Weight	25.7kDa(including fusion tag)
pI	6.9
Activity	NA
Storage	Storage buffer: 1XPBS. Store at -80°C and avoid repeated freeze-thaw cycles.
	kDa 40 30 25 20 15
	• <u>Interleukin-6</u> , a mental cytokine
Reference:	 Interleukin-6: from identification of the cytokine to development of targeted treatments Interleukin 6 in autoimmune and inflammatory diseases: a personal memoir Cardiovascular biology of interleukin-6 Tertiary structural models of human interleukin-6 and evaluation by comparison with X-ray and NMR structures The therapeutic potential of interleukin-6 in treating obesity Anti-interleukin-6 receptor antibody, tocilizumab, for the treatment of autoimmune diseases



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

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