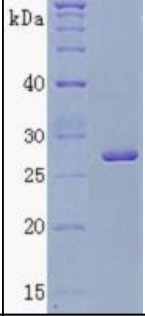


Recombinant Human IL6 Protein Data Sheet

Catalog #	hRP-A0301-EF011
Size	25 µg
Protein Name	Human interleukin 6
Protein Symbol	IL6
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
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.
	
Reference:	• Interleukin-6, a mental cytokine • 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™
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