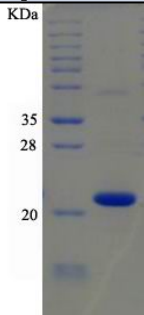


Recombinant Human IL1RN Protein Data Sheet

Catalog #	hRP-T2817-EF011
Size	25 µg
Protein Name	Human interleukin-1 receptor antagonist protein
Protein Symbol	IL1RN
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	BC009745.1
Uniprot Accession #	P18510-2
Description	Interleukin-1 receptor antagonist protein is a member of the interleukin 1 cytokine family. This protein inhibits the activities of interleukin 1, alpha (IL1A) and interleukin 1, beta (IL1B), and modulates a variety of interleukin 1 related immune and inflammatory responses.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	174aa(including fusion tag)
Molecular Weight	19.8kDa(including fusion tag)
pI	6.5
Activity	NA
Storage	Storage buffer: 20mM Tris.Cl, 50mM NaCl, 50% Glycerol, pH8.0. Store at -80°C and avoid repeated freeze-thaw cycles.
	
Reference:	• The efficacy and safety of interleukin-1-receptor antagonist anakinra in the treatment of systemic juvenile idiopathic arthritis • Schnitzler syndrome: response to anakinra in two cases and a review of the literature • The future of the IL-1 receptor antagonist anakinra: from rheumatoid arthritis to adult-onset Still's disease and systemic-onset juvenile idiopathic arthritis • Conformation and side chains environments of recombinant human interleukin-1 receptor antagonist (rh-IL-1ra) probed by raman, raman optical activity, and UV-resonance Raman spectroscopy • Interleukin 1 Receptor Antagonist Deficiency Presenting as Infantile Pustulosis Mimicking Infantile Pustular Psoriasis

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