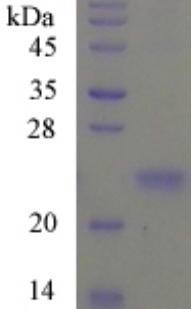


Recombinant Human CSF3 Protein Data Sheet

Catalog #	hRP-Z0547-EF041
Size	25 µg
Protein Name	Homo sapiens colony stimulating factor 3 (granulocyte)
Protein Symbol	CSF3
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	NM_000759.1
Uniprot Accession #	P09919
Description	Granulocyte/macrophage colony-stimulating factors are cytokines that act in hematopoiesis by controlling the production, differentiation, and function of 2 related white cell populations of the blood, the granulocytes and the monocytes-macrophages. This CSF induces granulocytes.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	none
Peptide Length	207aa
Molecular Weight	22.3kDa
pI	5.9
Activity	NA
Storage	Storage buffer: 20mM Tris.Cl, 50mM NaCl, 50% Glycerol, pH8.0. Store at -80°C and avoid repeated freeze-thaw cycles.
	



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