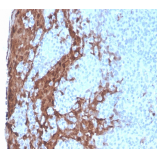


Granulocyte Colony Stimulating Factor Recombinant Monoclonal Mouse Antibody (rCSF3/900)



Product Description

This MAbs recognizes granulocyte-colony stimulating factor (G-CSF) in the cytoplasm of mature granulocytes. It shows no reactivity with any other cell types. Markers of myeloid cells are useful in the identification of different levels of cellular differentiation. It reacts with early precursor and mature forms of myeloid cells. It is useful for the detection of myeloid leukemias and granulocytic sarcomas. It can be used as a marker of granulocytes in normal tissues or inflammatory processes. G-CSF is a pleiotropic cytokine that influences differentiation, proliferation and activation of the neutrophilic granulocyte lineage. The human G-CSF cDNA encodes a 207 amino acid precursor containing a 29 amino acid signal peptide that is proteolytically cleaved to form a 178 amino acid residue mature protein. Two G-CSF s, which are identical except for a three amino acid deletion in the amino-terminus of one form of the protein have been isolated from human cells. Murine and human G-CSF s share 73% sequence identity at the amino acid level.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 3638, Anti-Granulocyte Colony Stimulating Factor (rCSF3/900)

Product attributes

Antibody number	#3638
Antibody reactivity (target)	Granulocyte Colony Stimulating Factor
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rCSF3/900
Isotype	IgG1, kappa
Molecular weight	19 kDa
Synonyms	Pluripoietin; Filgrastim; Lenograstim; CSF3OS; G-CSF; chromosome 17 open reading frame 33 (C17orf33); colony stimulating factor 3 (granulocyte)
Human gene symbol	CSF3
Entrez gene ID	1440
SwissProt	P09919
Unigene	2233
Immunogen	Recombinant full-length human CSF3 protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human, Macaque Monkey
Positive control	Tonsil or lymph node (IHC).
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Granulocytes, Myeloid cells
Tumor expression	Leukemia/lymphoma, Sarcoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on August 24, 2025 at 09:37:42 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/granulocyte-colony-stimulating-factor-recombinant-monoclonal-mouse-antibody-rcsf3-900/>