IL-6 Monoclonal Rat Antibody (MP5-20F3)

Product Description

IL-6 is a potent lymphoid cell growth factor that stimulates the growth and survivability of certain B-cells and T-cells. IL-6 plays a role in host defense, acute phase reactions, immune responses, and hematopoiesis. IL-6 is expressed by T-cells, B-cells, monocytes, fibroblasts, hepatocytes, endothelial cells and keratinocytes.

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 0663, Anti-IL-6 (MP5-20F3)

Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	0663		
Antibody reactivity (target)	IL-6		
Antibody type	Primary		
Host species	Rat		
Clonality	Monoclonal		
Clone	MP5-20F3		
Isotype	lgG1, kappa		
Molecular weight	21 kDa		
Synonyms	B cell differentiation factor; B-cell stimulatory factor 2 (BSF-2); CTL differentiation factor (CDF); Cytotoxic T cell differentiation factor; Hepatocyte stimulating factor; HPGF; HSF; Hybridoma growth factor Interferon beta-2; Hybridoma plasmacytoma growth factor; Interferon beta-2 (IFNB2)		
Entrez gene ID	16193 (Mouse)		
SwissProt	P08505 (Mouse)		
Unigene	1019 (Mouse)		
Immunogen	Recombinant mouse IL-6 protein		
Antibody target cellular localization	Secreted (extracellular)		
Species reactivity	Mouse		
Expected antibody applications	Flow (intracellular) (published for clone), Functional studies (published for clone)		
Antibody application notes	Flow cytometry: 0.5-1 ug/million cells in 0.1 mL, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Does not react with human or rat; others not known, Neutralization Studies order Ab without azide, Optimal dilution for a specific application should be determined by user		
Positive control	Mouse heart		
Positive control Shipping condition	Mouse heart Room temperature		
Shipping condition	Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Shipping condition Storage Conditions	Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when		
Shipping condition Storage Conditions Shelf life	Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended		
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% aside, Purified: BSA-free: 1 mg/mL in PBS/0.05% aside, Purified: 0.2 mg/mL in PBS/0.05% aside,		
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% 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		

Email: techsupport@biotium.com

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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

- 1. J Immunol (1990) 145(12): 4185-4191. (functional studies)
- 2. J Immunol (2013) 190:703-711. (functional studies; Flow)

This datasheet was generated on October 31, 2025 at 12:37:01 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-il-6-mouse-mp5-20f3/