



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)**

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)

Product attributes	
Antibody number	#0663
Antibody reactivity (target)	IL-6
Antibody type	Primary
Host species	Rat
Clonality	Monoclonal
Clone	MP5-20F3
Isotype	IgG1, 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
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Cytokines, Immunology, Inflammation

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

Dye Features
CF®405S Features
CF®488A Features
CF®568 Features
CF®594 Features
CF®640R Features
CF®647 Features

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 LI-COR Bioscience.