CD71 / Transferrin Receptor Monoclonal Mouse Antibody (66IG10)



Product Description

This antibody recognizes a ~90-95 kDa protein which is identified as cell surface transferrin receptor (CD71), a disulfide-bonded homodimeric glycoprotein of 180-190 kDa. This MAb is highly specific to CD71 and shows no cross-reaction with other related proteins. Its epitope is localized in the extracellular domain of CD71. Ligand for transferrin receptor is the serum iron transport protein, transferrin. This receptor is broadly distributed in carcinomas, sarcomas, leukemias, and lymphomas. CD71/Transferrin receptor has been reported to be associated with cell proliferation in both normal and neoplastic tissues and useful in predicting clinical behavior or response to therapy in a number of malignancies including breast cancer.

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 0871, Anti-CD71 (66IG10)

Product attribute

Call us: 800-304-5357

Product attributes				
Antibody number	#0871			
Antibody reactivity (target)	CD71, Transferrin Receptor			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	66IG10			
Isotype	IgG1, kappa			
Molecular weight	85-95 kDa (monomer); 190 kDa (dimer)			
Synonyms	Mtvr-1, p90, TFR1, TFRC transferrin receptor (p90 CD71), TRFR			
Human gene symbol	TFRC			
Entrez gene ID	7037			
SwissProt	P02786			
Unigene	529618			
Immunogen	Human thymocytes (T-cells)			
Antibody target cellular localization	Plasma membrane			
Verified antibody applications	Flow (surface) (verified)			
	Human			
Species reactivity	Human			
Species reactivity Expected antibody applications	Human Flow, surface (published for clone), IF (published for clone), IP (published for clone)			
Expected antibody	Flow, surface (published for clone), IF (published for clone), IP			
Expected antibody applications	Flow, surface (published for clone), IF (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Flow Cytometry 0.5-1 ug/mllilion cells/0.1 mL, Optimal dilution for a			
Expected antibody applications Antibody application notes	Flow, surface (published for clone), IF (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/ml., Flow Cytometry 0.5-1 ug/million cells/0.1 ml., Optimal dilution for a specific application should be determined by user CCRF-CEM, Jurkat or K-562. Human placenta or breast			
Expected antibody applications Antibody application notes Positive control	Flow, surface (published for clone), IF (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user CCRF-CEM, Jurkat or K-562. Human placenta or breast carcinoma.			
Expected antibody applications Antibody application notes Positive control Shipping condition	Flow, surface (published for clone), IF (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence 1-2 ug/ml. Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user CCRF-CEM, Jurkat or K-562. Human placenta or breast carcinoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Expected antibody applications Antibody application notes Positive control Shipping condition Storage Conditions	Flow, surface (published for clone), IF (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user CCRF-CEM, Jurkat or K-562. Human placenta or breast carcinoma. 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			
Expected antibody applications Antibody application notes Positive control Shipping condition Storage Conditions Shelf life	Flow, surface (published for clone), IF (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence 1-2 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user CCRF-CEM, Jurkat or K-562. Human placenta or breast carcinoma. 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			
Expected antibody applications Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	Flow, surface (published for clone), IF (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user CCRF-CEM, Jurkat or K-562. Human placenta or breast carcinoma. 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, Purffied: 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA/0.95% azide, Purffied; 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA/0.95% aride, Purffied; 0.85A-free: 1 mg/mL in PBS/0.05% BSA/0.05% azide, Purffied; 0.25 mg/mL in PBS/0.05%			
Expected antibody applications Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Flow, surface (published for clone), IF (published for clone), IP (published for clone) Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user CCRF-CEM, Jurkat or K-562. Human placenta or breast carcinoma. 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/0.55% azide, Purified, BSA-free: 1 mg/mL in PBS without azide			

Chinese hamster ovary cells. Inquire for the specific lot.

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,	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 (1999) 162: 1488-1495. (cell internalization; IF)
- 2. J Virol (2001) 75(24): 12347-12358. (Flow, surface)
- 3. PNAS ÙSA (2003) 100(20): 11583-11588. (IP)
- 4. Mol Biol Cell (2003) 14(12): 4745-5124. (cell internalization; IP)