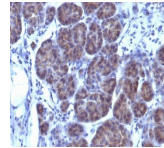


GLG1 Monoclonal Mouse Antibody (GLG1/970)



Product Description

This MAb recognizes a protein of 134 kDa, which binds fibroblast growth factor and E-selectin (cell-adhesion lectin on endothelial cells mediating the binding of neutrophils). Fucosylation is essential for binding to E-selectin. It contains sialic acid residues and 16 Cys-rich GLG1 repeats. This MAb can be used to stain the Golgi complex in cell or tissue preparations and can be used as a Golgi marker in subcellular fractions. It produces a diffuse staining pattern of the Golgi zone in normal and malignant cells. This MAb is an excellent marker for human cells in xenographic model research. It reacts specifically with human cells. The Golgi apparatus is an organelle present in all eukaryotic cells that forms a part of the endomembrane system. The primary function of the Golgi apparatus is to process and package macromolecules synthesized by the cell for exocytosis or use within the cell. The Golgi is made up of a stack of flattened, membrane-bound sacs known as cisternae, with three functional regions: the cis face, medial region and trans face. Each region consists of various enzymes that selectively modify the macromolecules passing through them, depending on where they are destined to reside. Several spherical vesicles that have budded off of the Golgi are present surrounding the main cisternae. The Golgi tends to be more pronounced and numerous in cells that make and secrete many substances such as plasma B cells. 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.

Product attributes

Antibody number	#0970
Antibody reactivity (target)	GLG1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	GLG1/970
Isotype	IgG1, kappa
Molecular weight	134 kDa
Synonyms	CFR-1; Cysteine-rich fibroblast growth factor receptor; E-selectin ligand-1 (ESL-1); Golgi Glycoprotein 1 (GLG1); Golgi apparatus protein 1; Golgi sialoglycoprotein MG-160; Slectin, endothelial cell, ligand
Human gene symbol	GLG1
Entrez gene ID	2734
SwissProt	Q92896
Unigene	109731
Immunogen	Golgi fraction from human liver cells
Antibody target cellular localization	Golgi apparatus
Verified antibody applications	IHC (FPPE) (verified)
Species reactivity	Human
Antibody application notes	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 mouse or rat, others not known. Immunohistology formalin-fixed 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes. Immunocytochemistry Acetone or paraformaldehyde fixed 0.5-1 ug/mL. Western blotting 0.5-1.0 ug/mL. Flow Cytometry 0.5-1.0 ug/million cells in 0.1 mL. Optimal dilution for a specific application should be determined by user
Positive control	HepG2, A431 or HeLa cells. Placenta, Tonsil, Testis and Ovary.
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	Organelle markers

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, LI-COR Bioscience.