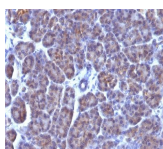


Golgi Complex Monoclonal Mouse Antibody (AE-6)



Product Description

This MAb recognizes an antigen associated with the Golgi complex in human cells only. It 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.

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 0868, Anti-Golgi Complex (AE-6)

Product attributes

Antibody number	#0868
Antibody reactivity (target)	Golgi Complex
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	AE-6
Isotype	IgG1, kappa
Molecular weight	Not Known
Synonyms	Golgi Complex; Golgi apparatus; Golgi zone
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	SU-DHL-1 large cell lymphoma cells.
Antibody target cellular localization	Golgi apparatus
Verified antibody applications	IHC (FFPE) (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
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.

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 November 22, 2024 at 10:27:01 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-golgi-complex-ae-6/>