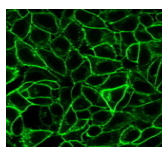


Endoglin / CD105 Monoclonal Mouse Antibody (ENG/1621)



Product Description

CD105/Endoglin is a Type I membrane glycoprotein located on cell surfaces and is part of the TGF-beta receptor complex. This protein has been found on endothelial cells, activated macrophages, fibroblasts, and smooth-muscle cells. Endoglin has a role in the development of the cardiovascular system and in vascular remodeling. Its expression is regulated during heart development. CD105 is highly expressed in endothelial cells during tumor angiogenesis and inflammation, with weak or negative expression in vascular endothelium of normal tissues. Angiogenesis is a promising prognostic marker in a variety of tumors. Endoglin is a more specific and sensitive marker for tumor angiogenesis than CD31 or CD34, as it labels only newly-formed blood vessels and may serve as a prognostic marker for prostate adenocarcinoma, and cancers of the lung, stomach, breast, and brain.

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 1621, Anti-CD105|Endoglin (ENG/1621)

Product attributes

Antibody number	#1621
Antibody reactivity (target)	CD105, Endoglin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ENG/1621
Isotype	IgG1, kappa
Molecular weight	84 kDa
Synonyms	CD105; END; Endoglin; Eng; HHT1; S-endoglin
Human gene symbol	ENG
Entrez gene ID	2022
SwissProt	P17813
Unigene	76753
Immunogen	Recombinant extracellular fragment of human Endoglin protein (exact sequence is proprietary)
Antibody target cellular localization	Plasma membrane
Species reactivity	Horse, Human, Monkey, Pig, Rat
Verified antibody applications	Flow (verified), IF (verified)
Antibody application notes	Flow cytometry: 1-2 ug/million cells, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 2-4 ug/mL, ELISA: For coating order antibody without BSA, For functional studies, order Ab without azide, Optimal dilution for a specific application should be determined by user
Positive control	HeLa, A431 and HL-60 cells. Colon.
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 research areas	Cancer
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
Cell/tissue expression	Endothelial cells
Tumor expression	Brain cancer, Breast cancer, Gastrointestinal cancer, Lung cancer, Prostate cancer
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 September 11, 2025 at 01:35:04 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/endoglin-cd105-monoclonal-mouse-antibody-eng-1621/>