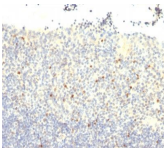


FOXP3 Monoclonal Mouse Antibody (3G3)



Product Description

Recognizes a protein of 47-55 kDa, which is identified as FOXP3. Its precise epitope is not known, but it has been mapped to the N-terminal portion (amino acids 2-136) of the protein. The FOX family of transcription factors is a large group of proteins that share a common DNA binding domain termed a winged-helix or forkhead domain. During early development, FOXP1 and FOXP2 are expressed abundantly in the lung, with lower levels of expression in neural, intestinal and cardiovascular tissues, where they act as transcription repressors. FOXP1 is widely expressed in adult tissues, while neoplastic cells often exhibit a dramatic change in expression level or localization of FOXP1. Mutations in FOXP3 gene cause IPEX, a fatal, X-linked inherited disorder characterized by immune dysregulation. The FOXP3 protein is essential for normal immune homeostasis. Specifically, FOXP3 represses transcription through a DNA binding forkhead domain, thereby regulating T cell activation. 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 0645, Anti-FOXP3 (3G3)**

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

- PNAS USA (2006) 103(17): 6659-6664. (Flow; epitope mapping)
- Cytometry Part A (2009) 75A: 1040-1050. (Flow; comparison of anti-FOXP3 clones and fixation methods)

Product attributes

Antibody number	#0645
Antibody reactivity (target)	FOXP3
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	3G3
Isotype	IgG1, kappa
Molecular weight	47-55 kDa
Synonyms	JM2, AIID, DIETER, IPEX, MGC141961, MGC141963, PIDX, XPID, FOXP3delta7, forkhead box protein P3, immune dysregulation, polyendocrinopathy, enteropathy, X-linked, immunodeficiency, polyendocrinopathy, enteropathy, X-linked, scurfir
Human gene symbol	FOXP3
Entrez gene ID	50943
SwissProt	Q9BZS1
Unigene	247700
Immunogen	Full-length His-tagged mouse FOXP3 protein
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Human, Monkey, Mouse
Expected antibody applications	Flow (intracellular) (published for clone)
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, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
Positive control	Tonsil, lymph node or Breast carcinoma
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	Cancer, Developmental biology, Endocrinology, Transcription factors

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.