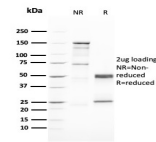


GATA-3 Monoclonal Mouse Antibody (GATA3/2446)



Product Description

GATA-3 is a zinc finger transcription factor and plays an important role in promoting and directing cell proliferation, development, and differentiation in many tissues and cell types. GATA-3 expression is primarily seen in breast and urothelial carcinomas. Therefore, GATA3 antibody can be useful in the identification of unknown primary carcinoma when carcinomas of the breast or bladder are a possibility. 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	#2446
Antibody reactivity (target)	GATA-3
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	GATA3/2446
Isotype	IgG2b, kappa
Molecular weight	~50 kDa
Synonyms	GATA3; GATA binding protein-3; GATA-binding factor 3; GATA3; HDR; HDRS; Transacting T-cell-specific transcription factor GATA-3
Human gene symbol	GATA3
Entrez gene ID	2625
SwissProt	P23771
Unigene	524134
Immunogen	Recombinant fragment of human GATA3 protein (around aa 357-436) (exact sequence is proprietary)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	MCF-7 or T47D cells. Breast or Bladder 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
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended

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
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, LI-COR Bioscience.

This datasheet was generated on July 23, 2024 at 10:31:58 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/gata-3-monoclonal-mouse-antibody-gata3-2446/>