



Glowing products for science

Trefoil factor 2 / TFF2 Monoclonal Mouse Antibody (GE16C)



Product Description

Trefoil peptides are protease resistant molecules secreted throughout the gut that play a role in mucosal healing. These peptides contain three intrachain disulfide bonds, forming the trefoil motif, or P-domain.

This antibody is available purified, with BSA and azide (0.2 mg/mL) or purified, BSA- and azide-free (1 mg/mL). **Catalog number key for antibody number 3060, Anti-TFF2|Trefoil factor 2 (GE16C)**

Call us : [800-304-5357](tel:800-304-5357) Email: techsupport@biotium.com

Product attributes

Antibody number	#3060
Antibody reactivity (target)	TFF2, Trefoil factor 2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	GE16C
Isotype	IgM, kappa
Molecular weight	14 kDa
Synonyms	SML1; SP; Spasmolysin; Spasmolytic polypeptide; TFF2; trefoil factor 2 (spasmolytic protein 1); Trefoil factor 2; Trefoil factor 2 precursor
Human gene symbol	TFF2
Entrez gene ID	7032
SwissProt	Q03403
Unigene	2979
Immunogen	Synthetic peptide corresponding to a 16-amino acid C-terminal sequence of the human Trefoil factor 2.
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Expected antibody applications	IHC (FFPE) (published for clone)
Antibody application notes	ELISA: For coating, order Ab without BSA & azide; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	Human stomach or duodenum tissues (IHC).
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
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
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.

References

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

Clin Cancer Res (2005) 11(18): 6472-6478. (IHC, FFPE)

This datasheet was generated on July 16, 2025 at 12:04:12 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/trefoil-factor-2-tff2-monoclonal-mouse-antibody-ge16c/>