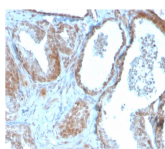


TIGIT / VSTM3 / VSIG9 Monoclonal Mouse Antibody (TIGIT/3106)



Product Description

TIGIT is a checkpoint inhibitor which binds with high affinity to the poliovirus receptor (PVR), causing increased IL10 secretion, decreased IL12B secretion. TIGIT binding to PVR also causes the suppression of T cell activation by promoting the generation of mature immuno-regulatory dendritic cells. It is expressed at low levels on natural killer (NK) cells, as well as peripheral memory and regulatory CD4 T cells. At the protein level, it is upregulated following the activation of these cells. Functionally, TIGIT is similar to CTLA4. The ligands for TIGIT include CD155 (signal abrogation) and CD226 (signal stimulation). It has been demonstrated to be upregulated on T cells in many cancers and is a immuno-oncology target for therapy.

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 3106, Anti-TIGIT|VSIG9|VSTM3 (TIGIT/3106)

Product attributes

Antibody number	#3106
Antibody reactivity (target)	TIGIT, VSIG9, VSTM3
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	TIGIT/3106
Isotype	IgG2b, kappa
Molecular weight of antigen	26 kDa
Synonyms	T-cell immunoreceptor with Ig and ITIM domains; V-set and immunoglobulin domain containing 9 (VSIG9); V-set and transmembrane domain containing 3 (VSTM3); Washington University Cell Adhesion Molecule (WUCAM)
Human gene symbol	TIGIT
Entrez gene ID	201633
SwissProt	Q4951
Unigene	421750
Immunogen	Recombinant fragment (around aa 22-141) of human TIGIT protein (exact sequence is proprietary)
Antibody target cellular localization	Plasma membrane
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Positive control	Tonsil or Prostate 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
Validated in protein array	Monospecific
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 (<i>Bos taurus</i>), 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 June 20, 2026 at 02:26:21 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/tigit-vstm3-vsigt9-mono-clonal-mouse-antibody-tigit-3106/>