

CD357 / GITR Monoclonal Rat Antibody (DTA-1)

Product Description

GITR (glucocorticoid-induced TNFR-related gene) is a member of the TNF-receptor superfamily, also known as TNFRSF18. It is expressed at low levels on resting T lymphocytes and at high levels on CD25 CD4 Tregs. The expression of GITR on T cells can be upregulated upon activation. Interaction of GITR with its ligand (GITRL) has been demonstrated to augment T cell activation, proliferation, cytokine production as well as MAPKs and NF- κ B activation, and abrogate the inhibitory function of CD25 CD4 Tregs. In vivo activation of GITR causes development of autoimmune diseases and restores the suppressed immune response. **Catalog number key for antibody number 0976, Anti-GITR|Tnfrsf18 (DTA-1)**

Product attributes

Antibody number	#0976
Antibody reactivity (target)	GITR, Tnfrsf18
Antibody type	Primary
Host species	Rat
Clonality	Monoclonal
Clone	DTA-1
Isotype	IgG2b, kappa
Molecular weight	66-70 kDa (Homodimer)
Synonyms	Activation-inducible TNF-related ligand; AITRL; GITRL; Glucocorticoid induced TNFR related protein ligand; Glucocorticoid-induced TNF-related ligand; hGITRL; TL6; TNFSF18; Tumor necrosis factor (ligand) superfamily member 18
Human gene symbol	TNFRSF18
Entrez gene ID	21936
SwissProt	O35714
Unigene	3180
Immunogen	Mouse CD25+ CD4+ T cells
Antibody target cellular localization	Plasma membrane
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone)
Species reactivity	Mouse
Antibody application notes	Flow cytometry: 0.5-1 ug/million cells in 0.1 mL, 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, Optimal dilution for a specific application should be determined by user
Positive control	CD25+CD4+Tregs.
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
Cell/tissue expression	T-cells
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 research areas	Immunology, Inflammation

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.

1. Clin Cancer Res (2013) 19(5): 1044-53. (functional studies)
2. J Immunol (2019) 203:864-872. (Flow)