CD90 / Thy1 Monoclonal Mouse Antibody (AF-9)

Product Description

Recognizes a protein of 18-35 kDa, identified as CD90 (also known as Thy1). CD90 is a member of the immunoglobulin superfamily. It may contribute to inhibition of proliferation/differentiation of hematopoietic stem cells and neuron memory formation in the CNS. It consists of a single Ig domain (112 amino acids; 25-35 kDa) inserted into the cell membrane via a GPI anchor. Expressed by hematopoietic stem cells and neurons in all species studied. Its highly expressed in connective tissue and various fibroblast and stromal cell lines, expressed on all thymocytes and peripheral T cells in mice, but in humans expressed only on small % fetal thymocytes, 10-40% of CD34 cells in bone marrow, and <1% of CD3 CD4 lymphocytes in peripheral circulation. It is also expressed by human lymph node HEV endothelium but not other endothelia. Lastly, it is expressed by a limited number of lymphoblastoid and leukemic cell lines.

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 0987, Anti-CD90 (AF-9)

Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	9987		
Antibody reactivity (target)	CD90, Thy1		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	AF-9		
Isotype	IgG1, kappa		
Molecular weight	18-35 kDa		
Synonyms	CD54; Cell surface glycoprotein P3.58; Human rhinovirus receptor; ICAM-1; Intercellular adhesion molecule 1; Ly 47; Major group rhinovirus receptor; MALA2; MyD10; Surface antigen of activated B cells		
Human gene symbol	THY1		
Entrez gene ID	7070		
SwissProt	P04216 644697		
Unigene	7070 P04216 644697 Human T-lymphoma cells Plasma membrane Cow, Human		
Immunogen	Human T-lymphoma cells		
Antibody target cellular localization	Plasma membrane		
Species reactivity	Cow, Human		
Expected antibody applications	Plasma membrane Cow, Human Flow, surface (published for clone), IHC (FFPE) (published for clone), IF (published for clone)		
Antibody application notes	Higher concentration may be required for direct detection using		
	clone), IF (published for clone) 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 MR-32, CCRF-CEM or MOLT-4 cells. Human uterus.		
Positive control	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mlL, Flow Cytometry 0.5-1 ug/mllion cells/0.1 ml_, Optimal dilution for a specific application should be determined by user IMR-32, CCRF-CEM or MOLT-4 cells. Human uterus.		
Positive control Shipping condition	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		
	secondary antibody, İmmunofluorescence: 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 IMR-32, CCRF-CEM or MOLT-4 cells. Human uterus.		
Shipping condition	secondary antibody, İmmunofluorescence: 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 IMR-32, CCRF-CEM or MOLT-4 cells. Human uterus. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Shipping condition Storage Conditions	secondary antibody, İmmunofluorescence: 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 IMR-32, CCRF-CEM or MOLT-4 cells. Human uterus. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when		
Shipping condition Storage Conditions Shelf life	secondary antibody, İmmunofluorescence: 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 IMR-32, CCRF-CEM or MOLT-4 cells. Human uterus. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended		
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	secondary antibody, İmmunofluorescence: 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 IMR-32, CCRF-CEM or MOLT-4 cells. Human uterus. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% 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, Purflect: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflect 0.2 mg/mL in PBS/0.05% azide, Purflect 0.2 mg/mL in PB		

Email: btinfo@biotium.com

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. Int J Stem Cell Res Ther (2018) 5:056. (IHC, FFPE)
- 2. Biology Open 2017 6: 1423-1433. (Flow)
- 3. Gastroenterol (2009) 136(2): P705-714. (IF)

This datasheet was generated on November 8, 2025 at 02:05:03 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-cd90-af-9/