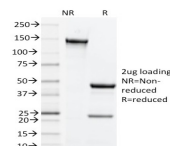


# CD134 (Mouse) Monoclonal Rat Antibody (OX-86)



## Product Description

CD134 is a type I integral membrane protein. This receptor is expressed on activated CD4 and CD8 T cells and B cells. The CD134 binds to CD134 ligand (CD252) to provide a costimulatory signal that is independent of CD28. CD134 is involved in coordinating CD4 T cell selection, migration and cytokine differentiation in T helper (Th)1 and Th2 cells. CD134 is also involved in the stimulation of T cells, T dependent humoral response and generation of optimal CD4 T cell responses in vivo and in vitro. CD134 is expressed on activated CD4 T lymphocytes, and its ligand, CD134L, is found preferentially on activated B cells. Engagement of CD134 with its ligand, CD134L, delivers a strong costimulatory signal to effector T cells.

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 2004, Anti-CD134, (OX-86)**

## Product attributes

Antibody number	#2004
Antibody reactivity (target)	CD134,
Antibody type	Primary
Host species	Rat
Clonality	Monoclonal
Clone	OX-86
Isotype	IgG1, kappa
Molecular weight	~50 kDa
Synonyms	CD134; IMD16; Lymphoid activation antigen (ACT35); OX40 cell surface antigen; OX40L receptor; TAX transcriptionally activated glycoprotein 1 receptor; Tumor necrosis factor receptor superfamily member 4 (TNFRSF4); Txgp11
Entrez gene ID	22163
SwissProt	P47741
Unigene	13885
Immunogen	Purified recombinant Mouse OX-40 antigen
Antibody target cellular localization	Plasma membrane
Species reactivity	Mouse
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone), IF (published for clone), WB (published for clone)
Positive control	MPBMC's. Spleen, Thymus or Lymph Node.
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.
Cell/tissue expression	B-cells, T-cells

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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. J Immunol (2000) 164:2160-2169. (functional studies)
2. Exp Med (2003) 198(5): 737-746. (Flow, surface)
3. Am J Respir Crit Care Med (2010) 181: 688-698. (WB)

J Leukoc Biol (2007) 81: 766-774. (IF)

This datasheet was generated on January 7, 2026 at 03:49:28 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/cd134-mouse-monoclonal-rat-antibody-ox-86/>