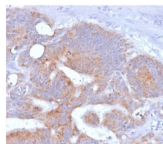


CD86 Recombinant Monoclonal Mouse Antibody (rC86/1146)



Product Description

This antibody recognizes a protein of 70 kDa, which is identified as CD86. CD86 is a type I transmembrane glycoprotein and a member of the immunoglobulin superfamily of cell surface receptors. It is expressed at high levels on resting peripheral monocytes and dendritic cells and at very low density on resting B and T lymphocytes. CD86 expression is rapidly upregulated by B cell specific stimuli with peak expression at 18 to 42 hours after stimulation. CD86, along with CD80/B71, is an important accessory molecule in T cell co-stimulation via its interaction with CD28 and CD152/CTLA4. Since CD86 has rapid kinetics of induction, it is believed to be the major CD28 ligand expressed early in the immune response. It is also found on malignant Hodgkin and Reed Sternberg (HRS) cells in Hodgkin's disease.

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 2083, Anti-CD86 (rC86/1146)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#2083
Antibody reactivity (target)	CD86
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rC86/1146
Isotype	IgG1, kappa
Molecular weight	70 kDa
Synonyms	Activation B7-2 antigen; B lymphocyte activation antigen B72; CD28 antigen ligand 2; CD28LG2; CLS1; CTLA-4 counter-receptor B7.2; Early T-cell co-stimulatory molecule 1; ETC-1; FUN-1; LAB72; Ly-58; MB7-2; TS/A-2
Human gene symbol	CD86
Entrez gene ID	942
SwissProt	P42081
Unigene	171182
Immunogen	Recombinant full-length human CD86 protein
Antibody target cellular localization	Plasma membrane
Verified antibody applications	Flow (verified), IF (verified), IHC (FFPE) (verified)
Species reactivity	Human, Mouse, Rat
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunohistology (formalin): 2-4 ug/mL for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris buffer with 1 mM EDTA pH 9.0 for 10-20 minutes followed by cooling at RT for 20 minutes. Optimal dilution for a specific application should be determined by user
Positive control	293T or Jukat cells. Monocytes and dendritic cells, activated T, B and natural killer cells in lymph node or tonsil.
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Cancer immunotherapy/immune checkpoint, Immunology
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	Dendritic cells, Monocytes/macrophages

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 January 28, 2026 at 12:54:45 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cd86-recombinant-monoclonal-mouse-antibody-rc86-1146/>