

# CD26 Monoclonal Mouse Antibody (202-36)

## Product Description

Recognizes a glycoprotein of 110 kDa, identified as CD26 (Workshop VI; Code: N-L039). It is an atypical serine protease belonging to the prolyl oligopeptidase family. It is expressed on lymphocyte cells and is upregulated during T-cell activation. CD26 is also expressed on activated B cells and natural killer cells and abundantly on epithelia. CD26 is implicated in a variety of biological functions including T-cell activation, cell adhesion with extracellular matrix such as fibronectin or collagens, and in HIV infection. Cross-linking of CD26 using this antibody dramatically enhances the anti-CD3-induced IL-2 production. In Western blotting, this MAb reacts with only glycosylated CD26, but not with the deglycosylated form. It does not prevent ADA binding to CD26.

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 0888, Anti-CD26 (202-36)**

## Product attributes

Antibody number	#0888
Antibody reactivity (target)	CD26
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	202-36
Isotype	IgG2b, kappa
Molecular weight	~110 kDa
Synonyms	ADA-binding protein (ADABP); Adenosine deaminase complexing protein 2 (ADCP-2); Dipeptidyl peptidase 4; Dipeptidyl peptidase 4 soluble form; Dipeptidyl peptidase IV membrane form; Dipeptidyl peptidase IV soluble form; Dipeptidyl peptidase, intestinal; Dipeptidylpeptidase 4; Dipeptidylpeptidase IV; DPP4; DPPIV; Intestinal dipeptidyl peptidase; T-cell activation antigen CD26; TP103
Human gene symbol	DPP4
Entrez gene ID	1803
SwissProt	P27487
Unigene	368912
Immunogen	Human T cell clone
Antibody target cellular localization	Plasma membrane
Species reactivity	Human, Rat
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone), IF (published for clone)
Antibody application notes	Immunohistology Frozen Only 0.5-1.0 ug/mL, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Does not react with pig or sheep, others not tested. Optimal dilution for a specific application should be determined by user
Positive control	YT, HEP-G2 cells, and lymphocytes. Lymph nodes and tonsils.
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	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	Epithelial 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. Cancer Res (2005) 65(15): 6950-6. (functional studies)
2. J Biol Chem (2001) 276(22): 19532-19539. (Flow)
3. BMC Cancer (2011) 11:51. (IF)

This datasheet was generated on January 28, 2026 at 02:26:34 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cd26-202-36/>