

CD37 Monoclonal Mouse Antibody (IPO-24)

Product Description

Recognizes a protein of 33-55 kDa, identified as CD37 (Workshop V; Code CD37.7). CD37 is strongly expressed on normal and neoplastic mature (slg) B-lymphocytes. In B-cell ontogeny, CD37 appears after the pre-B-cell stage, is maintained during peripheral B-cell development and is lost upon terminal differentiation into plasma cells.¹ CD37 is also present, at low densities, on resting and activated T cells, neutrophils, monocytes, and some myelomonocytic leukemia cells. It is absent from platelets, erythrocytes. CD37 is a member of a family of tetraspan transmembrane proteins, including CD9, CD53, CD63, CD81, and CD82. It associates other tetraspan transmembrane proteins and MHC class II molecules to form a large complex at the surface of B cells and play a role in signal transduction. CD37 is a valuable and stable marker for peripheral mature B-cells and corresponding malignancies like B-cell chronic lymphocytic leukemia (B-CLL), hairy cell leukemia (HCL), and all types of B-cell non-Hodgkin" lymphoma (B-NHL).

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 0170, Anti-CD37 (IPO-24)

Product attributes

Antibody number	#0170
Antibody reactivity (target)	CD37
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	IPO-24
Isotype	IgG2b, kappa
Molecular weight	33-55 kDa
Synonyms	Cell differentiation antigen 37; GP52 40; Tetraspanin-26; Tspan-26; TSPAN26
Human gene symbol	CD37
Entrez gene ID	951
SwissProt	P11049
Unigene	166556
Immunogen	Spleen cells of a patient with hairy cell leukemia
Antibody target cellular localization	Plasma membrane
Verified antibody applications	Flow (verified)
Species reactivity	Human
Expected antibody applications	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 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
Positive control	Daudi, Raji, Namalva, Ramos cells and human lymphocytes. Human tonsil and lymph node.
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
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	Lymphocytes, Myeloid cells
Tumor expression	Leukemia/lymphoma

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.

- Blood (2010) 116(7): 1025-1034. (IHC, FFPE; IF)
- PLoS One 12(10): e0185940. (Flow, surface; IF)