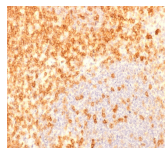


CD27 Recombinant Monoclonal Mouse Antibody (rLPFS2/1611)



Product Description

This antibody recognizes a protein of a disulfide-linked 120 kDa dimer, identified as CD27. It is expressed on the majority of peripheral T cells, medullary thymocytes, memory-type B cells, and natural killer cells. It is a transmembrane phosphoglycoprotein that belongs to the tumor necrosis factor receptor (TNFR) superfamily. CD27 binds to its ligand CD70, a member of the TNF family, and induces T-cell co-stimulation and B-cell activation. It also interacts with TRAFs and is involved in activation of NFκB and SAPK/JNK and induces apoptosis. 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 1976, Anti-CD27 (rLPFS2/1611)**

Product attributes

Antibody number	#1976
Antibody reactivity (target)	CD27
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rLPFS2/1611
Isotype	IgG1, kappa
Molecular weight	120 kDa
Synonyms	LPFS2; S152; T cell activation antigen S152; T-cell activation antigen CD27; T14; TNFRSF7; TNFSF7; Tp55; Tumor necrosis factor receptor superfamily member 7
Human gene symbol	CD27
Entrez gene ID	939
SwissProt	P26842
Unigene	355307
Immunogen	Recombinant human CD22 protein (exact sequence is proprietary)
Verified antibody applications	Flow (verified), IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	Ramos cells. Human peripheral blood 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	Immunology
Cell/tissue expression	Lymphocytes

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
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, LI-COR Bioscience.