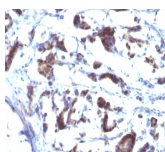


ZFYVE28 Monoclonal Mouse Antibody (LST2/2426)



Product Description

ZFYVE28 belongs to the Ist-2 family. It contains 1 FYVE-type zinc finger. The FYVE-type zinc finger mediates the interaction with phosphatidylinositol 3-phosphate (PI3P) and localization to early endosome membranes when not mono-ubiquitinated at Lys-87. Mono-ubiquitination at Lys-87 prevents binding to phosphatidylinositol 3-phosphate (PI3P) and localization to early endosome membranes. ZFYVE28 is a negative regulator of epidermal growth factor receptor (EGFR) signaling. It acts by promoting EGFR degradation in endosomes when not mono-ubiquitinated. The FYVE domain has been identified in a number of unrelated signaling molecules. This protein functions to recruit SMAD2 to the transforming growth factor-beta receptor. The FYVE domain is required to maintain the normal localization of this protein but is not involved in mediating interaction with SMADs.

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 2426, Anti-ZFYVE28 (LST2/2426)

Product attributes

Antibody number	#2426
Antibody reactivity (target)	ZFYVE28
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	LST2/2426
Isotype	IgG2a, kappa
Molecular weight	96 kDa
Synonyms	Lateral signaling target protein 2 homolog; LST2; LYST2; Zinc finger FYVE domain-containing protein 28 (ZFYVE28)
Human gene symbol	ZFYVE28
Entrez gene ID	57732
SwissProt	Q9H9
Unigene	292056
Immunogen	Recombinant full-length human ZFYVE28 protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Cytoplasmic
Species reactivity	Human
Positive control	PC3 cells. Testis, Brain, Colon, Kidney or Stomach.
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
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
Validated in protein array	Monospecific
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.

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 July 13, 2025 at 09:45:44 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/zfyve28-monoclonal-mouse-antibody-lst2-2426/>