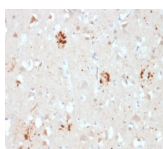


Ubiquitin Monoclonal Mouse Antibody (UBB/2122)



Product Description

Ubiquitin is a highly conserved and plays an essential role in the ubiquitin-proteasome pathway. In ubiquitination process, it is first activated by forming a thiol-ester complex with the activation component E1, which is then transferred to ubiquitin-carrier protein E2, followed by to ubiquitin ligase E3 for final delivery to epsilon-NH2 of the target protein lysine residue. IκB, p53, cdc25A, Bcl-2 etc. are shown as targets of ubiquitin-proteasome process as part of regulation of cell cycle progression, differentiation, cell stress response, and apoptosis. Moreover, ubiquitin have been reported to bind covalently with pathological inclusions which are resistant to degradation e.g. neurofibrillary tangles/paired helical filaments in Alzheimer's disease, Lewy bodies seen in Parkinson's disease, and Pick bodies found in Pick's disease etc.

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 2122, Anti-Ubiquitin (UBB/2122)

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

Product attributes

Antibody number	#2122
Antibody reactivity (target)	Ubiquitin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	UBB/2122
Isotype	IgG1, kappa
Molecular weight of antigen	9 kDa
Synonyms	Epididymis secretory protein LI 50; Polyubiquitin B; RPS27A; UBA52; UBA80; UBB; UBC; UBCEP1; UBCEP2; Ubiquitin; Ubiquitin B
Human gene symbol	UBB
Entrez gene ID	7314
SwissProt	P0CG47; P0CG48; P62979; P62987 and P62988
Unigene	356190 and 520348
Immunogen	Recombinant fragment of human Ubiquitin protein (around aa 1-119) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Cytoplasmic, Plasma membrane, Nucleus
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, ELISA For coating, order Ab without BSA. Optimal dilution for a specific application should be determined by user
Positive control	HeLa or Jurkat cells, Alzheimer s Brain
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
Validated in protein array	Monospecific
Antibody research areas	Ubiquitination
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 June 26, 2026 at 04:50:25 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/ubiquitin-monooclonal-mouse-antibody-ubb-2122/>