## **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. IkB, 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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2122, Anti-Ubiquitin (UBB/2122)

## Product attributes

| 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   | 9 kDa  |  |  |  |
| Synonyms   | Epididymis secretory protein Li 50; Polyubiquitin B; RPS27A;<br>UBA52; UBA80; UBB; UBC; UBCEP1; UBCEP2; Ubiquitin;<br>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<br>localization   | Cytoplasmic, Plasma membrane, Nucleus  |  |  |  |
|  | Human  |  |  |  |
| Species reactivity   | Human  |  |  |  |
| Species reactivity  Antibody application notes   | Human  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  |  |  |  |
|  | 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   |  |  |  |
| 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  |  |  |  |
| Antibody application notes  Positive control   | 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   |  |  |  |
| Antibody application notes  Positive control Shipping condition  | 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  HeLa or Jurkat cells. Alzheimer s Brain  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,   |  |  |  |
| Antibody application notes  Positive control Shipping condition Storage Conditions   | 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  HeLa or Jurkat cells. Alzheimer s Brain  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when  |  |  |  |
| Antibody application notes  Positive control Shipping condition Storage Conditions Shelf life  | 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  HeLa or Jurkat cells. Alzheimer s Brain  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  |  |  |  |
| Antibody application notes  Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate   | 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  HeLa or Jurkat cells. Alzheimer s Brain  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  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, BSA-free: 1 mg/mL in PBS/0.05% BSA, Purified; D.2 mg/mL in PBS/0.05% BSA, Purified; BSA-free: 1 mg/mL in PBS/0.05% BSA, Purified; BSA-free: 1 mg/mL in PBS/0.05% BSA/0.05% azide, Purified; D.2 mg/mL in PBS/0.05% BSA/0.05% azid |  |  |  |
| Antibody application notes  Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation                             | 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  HeLa or Jurkat cells. Alzheimer s Brain  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  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, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified; BSA-free: 1 mg/mL in PBS without azide  |  |  |  |
| Antibody application notes  Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation  Validated in protein array | 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  HeLa or Jurkat cells. Alzheimer s Brain  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  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 without azide  Monospecific  |  |  |  |

Call us: 800-304-5357 Email: btinfo@biotium.com

| Antibody # prefix | Conjugation           | Ex/Em (nm) | Laser line | Detection channel           | Dye Features     |
|-------------------|-----------------------|------------|------------|-----------------------------|------------------|
| BNC04             | CF®405S               | 404/431    | 405        | DAPI (microscopy),<br>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,<br>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 November 3, 2025 at 04:49:21 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/ubiquitin-monoclonal-mouse-antibody-ubb-2122/">https://biotium.com/product/ubiquitin-monoclonal-mouse-antibody-ubb-2122/</a>