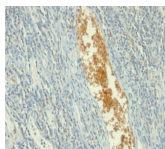


Adenosine Monophosphate Deaminase-3 Monoclonal Mouse Antibody (AMPD3/901)



Product Description

Adenosine Monophosphate Deaminase, isoform E (AMPD3) is a highly regulated enzyme that catalyzes the hydrolytic deamination of adenosine monophosphate to inosine monophosphate, a branch point in the adenylate catabolic pathway. AMPD3 gene encodes the erythrocyte (E) isoforms, whereas other family members encode isoforms that predominate in muscle (M) and liver (L) cells. It has 767 amino acids, a MW of ~ 90 kDa, and is assigned an EC 3.5.4.6. This MAb shows reactivity with cells of the erythroid lineage at all stages of maturation in the peripheral blood, bone marrow, and fetal liver. Non-erythroid lineages are negative by flow cytometry. This MAb is useful in the diagnosis of erythroleukemia, identification of bone marrow erythroid precursors, gating erythroid nucleated precursor cells from malignant cells in bone marrow specimens.

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 0901, Anti-Adenosine Monophosphate Deaminase-3 (AMPD3/901)

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

Product attributes

| | |
|---------------------------------------|--|
| research-areas | Cancer, Metabolism |
| Antibody number | #0901 |
| Antibody reactivity (target) | Adenosine Monophosphate Deaminase-3 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | AMPD3/901 |
| Isotype | IgG2b, kappa |
| Molecular weight | 90 kDa |
| Synonyms | AMPD3; Erythrocyte AMP Deaminase; AMP Deaminase 3 |
| Human gene symbol | AMPD3 |
| Entrez gene ID | 272 |
| SwissProt | Q01432 |
| Unigene | 501890 |
| Immunogen | Recombinant full-length human AMPD3 protein |
| Verified antibody applications | IHC (FFPE) (verified) |
| Antibody target cellular localization | Plasma membrane, Nucleus |
| Species reactivity | Human |
| Antibody application notes | Immunohistochemistry formalin-fixed 2-4 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Immunocytochemistry 1-2.0 ug/mL, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 1-2 ug/million cells in 0.1 mL, Optimal dilution for a specific application should be determined by user |
| Positive control | RBC. Fetal liver; Spleen or Placenta |
| 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, Metabolism |
| 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 | Red blood cells |
| Tumor expression | Erythroid cancer |

| 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 January 1, 2026 at 07:03:38 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-adenosine-monophosphate-deaminase-3-ampd3-ampd3901/>