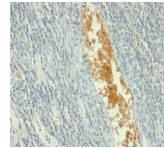


## Adenosine Monophosphate Deaminase-3 Monoclonal Mouse Antibody (AMPD3/901)



### 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](mailto: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)**

### 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	Immunohistology 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
Cell/tissue expression	Red blood cells
Tumor expression	Erythroid cancer

## Catalog Number Key for Primary Antibody Conjugates

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
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	

This datasheet was generated on January 30, 2023 at 07:33:50 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/monoclonal-anti-adenosine-monophosphate-deaminase-3-ampd3-ampd3901/>

