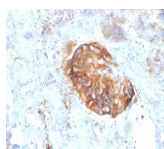


# MTAP Recombinant Monoclonal Mouse Antibody (rMTAP/1813)



## Product Description

Recognizes a protein of 31 kDa, which is identified as MTAP (5'-deoxy-5'-methylthioadenosine phosphorylase). It catalyzes the reversible phosphorolysis of methylthioadenosine, which is important in polyamine metabolism and for the salvage of adenine and methionine. The gene encoding MTAP is linked to the tumor suppressor gene, p16INK4A. Deficient levels of MTAP can occur in cancers primarily through co-deletion of the MTAP gene and the p16INK4A gene. Cells expressing MTAP and possessing adenine salvage pathway activity may be less susceptible to malignancy due to growth-inhibitory actions of agents (e. g. antifolates), whose mechanism of action, in part, involves this de novo purine pathway.

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 3646, Anti-MTAP (rMTAP/1813)**

## Product attributes

Antibody number	#3646
Antibody reactivity (target)	MTAP
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rMTAP/1813
Isotype	IgG1, kappa
Molecular weight	31 kDa
Synonyms	BDMF; DMSFH; DMSMFH; Epididymis luminal protein 249; HEL249; LGMBF; MeSAAdo phosphorylase; Methylthioadenosine phosphorylase; MSAP; MTA phosphorylase; MTAPase; S-methyl-5'-thioadenosine phosphorylase
Human gene symbol	MTAP
Entrez gene ID	4507
SwissProt	Q13126
Unigene	193268
Immunogen	Recombinant human MTAP protein fragment (aa97-196) (exact sequence is proprietary)
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Cytoplasmic, Nucleus
Species reactivity	Human
Positive control	HepG2, A431, HeLa or MCF-7 cells. Human kidney tissues (IHC).
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
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	<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>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 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	

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 April 3, 2025 at 08:37:58 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mtap-recombinant-monoclonal-mouse-antibody-rmtap-1813/>