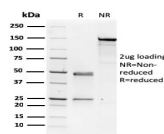


NME1 / nm23-H1 / NDPK-A Monoclonal Mouse Antibody (CPTC-NME1-2)



Product Description

The nm23 gene, a potential suppressor of metastasis, was originally identified by differential hybridization between two murine melanoma sub-lines, one with a high and the second with a low metastatic capacity. Highly metastatic sub-lines exhibit much lower levels of nm23 than less metastatic cells. Based on sequence analysis, nm23 appears highly related to nucleotide diphosphate kinases (NDP-K). In humans, NDP kinase A and B are identical to two isotypes of human nm23 homologs, namely nm23-H1 and H2, respectively. nm23-H2 is identical in sequence to PuF, a transcription factor that binds to nuclease hypersensitive elements at positions 142 to 115 of the human C-Myc promoter.

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 2247, Anti-NDPK-A|nm23-H1|NME1 (CPTC-NME1-2)

Product attributes

Antibody number	#2247
Antibody reactivity (target)	NDPK-A, nm23-H1, NME1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-NME1-2
Isotype	IgG2a, kappa
Molecular weight	23 kDa
Synonyms	AWD; GAAD; Granzyme A activated DNase; Metastasis inhibition factor NM23; NB; NBS; NDP kinase A; NM23 long variant; included; NM23H1B; NME/NM23 nucleoside diphosphate kinase 1; NME1-NME2 spliced read-through transcript; included; Non-metastatic protein 23; homolog 1; Nucleoside diphosphate kinase A; Tumor metastatic process-associated protein
Human gene symbol	NME1
Entrez gene ID	4830
SwissProt	P15531
Unigene	463456
Immunogen	Recombinant full-length human NME1 protein
Verified antibody applications	WB (verified)
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Human
Antibody application notes	Western blot 0.5-1.0 ug/mL; Optimal dilution for a specific application should be determined.
Positive control	HepG2 and HeLa whole cell lysates; heart, brain, placenta and lung carcinoma.
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 (<i>Bos taurus</i>), 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 November 27, 2024 at 07:39:10 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/nme1-nm23-h1-ndpk-a-mono-clonal-mouse-antibody-cptc-nme1-2/>