NME2 / nm23-H2 / NDPK-B Monoclonal Mouse Antibody (CPTC-NME2-2)



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). In humans, NDP kinases 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 promotor.

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 2248, Anti-NDPK-B|nm23-H2|NME2 (CPTC-NME2-2)

Product attributes

Product attributes			
Antibody number	#2248		
Antibody reactivity (target)	NDPK-B, nm23-H2, NME2		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	CPTC-NME2-2		
Isotype	IgG2a, kappa		
Molecular weight	17 kDa		
Synonyms	C myc purine binding transcription factor PUF; C-myc purine-binding transcription factor PUF; epididymis secretory sperm binding protein Li 155an; HEL-S-155an; Histldine protein kinase NDKB;; NDP kinase B; NDPKB; nm23-H2; NM23B; NME/NM23 nucleoside diphosphate kinase 2; nme2; Nucleoside diphosphate kinase B;		
Human gene symbol	NME2		
Entrez gene ID	4831		
SwissProt	P22392		
Unigene	463456		
Immunogen	Recombinant full-length human NME2 protein		
Verified antibody applications	IHC (FFPE) (verified), WB (verified)		
Antibody target cellular localization	Nucleus & cytoplasm		
Species reactivity	Human		
Positive control	PC3, A549, HeLa, Jurkat cells. Ubiquitously expressed in all tissues.		
Shipping condition	Room temperature		
Storage Conditions	Store at 2 to 8 $^{\circ}$ C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ 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		
Validated in protein array	Monospecific		
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.		

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), 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 5, 2025 at 02:17:55 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/nme2-nm23-h2-ndpk-b-monoclonal-mouse-antibody-cptc-nme2-2/