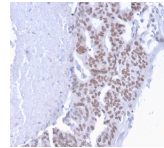


CHD4 Monoclonal Mouse Antibody (3F2/4)



Product Description

CHD4/Mi-2b is a component of the nucleosome remodeling and deacetylase (NuRD) complex which is a multi-subunit protein complex containing both histone deacetylase and nucleosome-dependent ATPase subunits. Current models predict that this complex function primarily in transcriptional repression. Accumulating evidence indicates that NuRD may regulate the transcription of specific genes by interacting with specific transcriptional factors. In addition, the NuRD complex may also participate in genome-wide transcriptional regulation through an association with histone tails. 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.

Product attributes

Antibody number	#3077
Antibody reactivity (target)	CHD4
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	3F2/4
Isotype	IgG2a, kappa
Molecular weight	238 kDa
Synonyms	ATP dependent helicase CHD4; Chromodomain helicase DNA binding protein 4; Mi 2 autoantigen 218 kDa protein; Mi 2b; Mi2 beta
Human gene symbol	CHD4
Entrez gene ID	1108
SwissProt	Q14839
Unigene	162233
Immunogen	Synthetic peptide (N- and C- terminal peptides were used for immunisation: 1. ASGLGSPSPCSAGSEEDM & 2. CSRLANRAPEPPPPQVAQQQ) – which peptide the antibody recognises has not been tested.
Verified antibody applications	IF (verified), IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Human, Mouse
Positive control	HeLa whole cell lysate or human placenta tissue lysate. Human breast 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

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
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, LI-COR Bioscience.