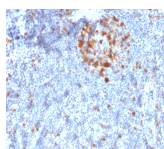


p34 / cdk1 Monoclonal Mouse Antibody (A17.1.1)



Product Description

Recognizes a 34 kDa protein (cdk1), identified as p34cdc2 (a catalytic subunit of Maturation Promoting Factor). Its epitope maps near the C-terminus of the protein and its core is thought to be LGTPNNEV (aa220-227 in murine cdc2). It shows no cross reaction with cdk2 p32. It supports the kinase activity. p34cdc2 plays a crucial role during cell division and is most active during mitosis. It is a serine/threonine kinase, which is activated by cyclin, by dephosphorylation of tyrosine residues. p34cdc2 is inactivated by a tyrosine kinase. This MAb reportedly inhibits the activation of p34cdc2 kinase by cyclins.

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 1083, Anti-Cdk1|p34cdc2 Serine-Threonine Kinase (A17.1.1)

Product attributes

Antibody number	#1083
Antibody reactivity (target)	Cdk1, p34cdc2 Serine-Threonine Kinase
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	A17.1.1
Isotype	IgG2a, kappa
Molecular weight	34 kDa
Synonyms	Cdc2; CDC28A; CDK1; CDKN1; CELL CYCLE CONTROLLER CDC2; Cell division control protein 2 homolog; Cell division cycle 2 G1 to S and G2 to M; Cell division protein kinase 1; Cyclin-dependent kinase 1; p34 Protein Kinase; P34CDC2
Human gene symbol	CDC2
Entrez gene ID	983
SwissProt	P06493
Unigene	334562
Immunogen	C-Terminal 2/3rds of Xenopus cdc2 expressed in E.coli
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Chicken, Guinea pig, Human, Mouse, Rat, Woodchuck, Xenopus laevis
Positive control	HeLa cells. Human tonsil (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 (<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 February 11, 2025 at 11:18:04 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/p34-cdk1-monooclonal-mouse-antibody-a17-1-1/>