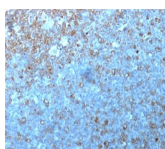


# p34 / cdk1 Monoclonal Mouse Antibody (POH-1)



## Product Description

Recognizes a 34 kDa protein, which is identified as cyclin dependent kinase 1 (cdk1) or p34cdc2 protein kinase. cdk1 / p34cdc2 plays a crucial role during cell division and is most active during mitosis. It is predominantly localized in the nucleus. It is a serine/threonine kinase, which is activated by cyclin, presumably by de-phosphorylation of tyrosine residues. Activated cdk1 / p34cdc2 performs specific functions during mitosis, including nuclear envelope breakdown and chromosome condensation.

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 0853, Anti-Cdk1 (POH-1)**

## Product attributes

Antibody number	#0853
Antibody reactivity (target)	Cdk1, p34
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	POH-1
Isotype	IgG1, 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	CDK1
Entrez gene ID	983
SwissProt	P06493
Unigene	334562
Immunogen	Purified recombinant human p34cdc2 fusion protein
Antibody target cellular localization	Nucleus & cytoplasm
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Species reactivity	Cow, Human, Mink, Monkey
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL. Does not react with mouse, rat, Xenopus, Drosophila, S. pombe, or S. cerevisiae; others not known. Immunohistology formalin-fixed 2-4 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user
Positive control	HeLa or MCF-7 Cells or Tonsil
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Cell cycle, Tumor suppressors
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 December 23, 2024 at 03:38:52 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-p34-cdk1-poh-1/>