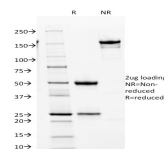


# Cdk2 / p34cdc2 Serine-Threonine Kinase Monoclonal Mouse Antibody (AN4.3)



## Product Description

In vertebrates, as in yeast, multiple cyclins have been identified, including a total of eight such regulatory proteins in mammals. In contrast to the situation in yeast, the Cdc2 p34 kinase is not the only catalytic subunit identified in vertebrates that can interact with cyclins. While Cdc2 p34 is essential for the G2 to M transition in vertebrate cells, a second Cdc2-related kinase has also been implicated in cell cycle control. This protein, designated cyclindependent kinase 2 (Cdk2), also binds to cyclins and its kinase activity is temporally regulated during the cell cycle. Several additional Cdc2-related cyclin dependent kinases have been identified. These include Cdk3, Cdk4, Cdk5, PCTAIRE-1, PCTAIRE-2, PCTAIRE-3, Cdk6 Cdk7, Cdk8 and KKIALLRE.

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 2316, Anti-Cdk2|p34cdc2 Serine-Threonine Kinase (AN4.3)**

## Product attributes

<b>Antibody number</b>	#2316
<b>Antibody reactivity (target)</b>	Cdk2, p34cdc2 Serine-Threonine Kinase
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	AN4.3
<b>Isotype</b>	IgG2a, kappa
<b>Molecular weight</b>	34 kDa
<b>Synonyms</b>	Cdc2 related protein kinase; CDC28; CDC2A; CDK1; CDK2; CDKN2; Cell division Kinase 2; Cell division protein kinase 2; cyclin dependent kinase 2 alpha; kinase Cdc2; MPF; p33 protein kinase
<b>Human gene symbol</b>	CDK2
<b>Entrez gene ID</b>	1017
<b>SwissProt</b>	P24941
<b>Unigene</b>	19192
<b>Immunogen</b>	Recombinant full-length human Cdk2 protein
<b>Antibody target cellular localization</b>	Nucleus
<b>Species reactivity</b>	Human, Mouse, Xenopus laevis
<b>Expected antibody applications</b>	WB (published for clone)
<b>Antibody application notes</b>	Inhibits activation of p34cdc2 kinase by cyclins; Immunoprecipitation; Kinase Assay; Western blot; Flow cytometry: 1-2 ug/mL; Optimal dilution for a specific application should be determined.
<b>Positive control</b>	HeLa cells, HeLa whole cell lysate.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	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.

## References

Note: References for this clone sold by other suppliers may be listed for expected applications.

Cell Cycle (2016) 15(23):3203-3209. (WB)

This datasheet was generated on November 27, 2024 at 12:30:34 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cdk2-p34cdc2-serine-threonine-kinase-monoclonal-mouse-antibody-an4-3/>