

# Cig2 Monoclonal Mouse Antibody (3A11/5)



## Product Description

Cig2 is a cyclin that functions as the catalytic subunit of cyclin-dependent kinases (Cdks) in the fission yeast *Schizosaccharomyces pombe*. It is a B-type, S phase cyclin that is required for both the G1/S and G2/M cell cycle transitions and is expressed in a sharp spike that peaks during the G1/S period. Cig2 binds to Cdc2 (Cdk), and the resulting Cdc2/Cig2 complex controls the G1/S transition of the cell cycle. Disruption of Cig2 delays the onset of mitosis. The expression of the Cig2 gene is dependent on Mlu1-binding factor (MBF), and the protein is destroyed during anaphase by the APC/cyclosome (APC/C) and Skp1/Cullin-1/F-box (SCF), thereby ensuring the spike expression pattern of Cig2. SCF regulates Cig2 levels in a dual manner, transcriptionally and post-translationally, while APC/C only destroys the protein.

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 2827, Anti-Cig2 (3A11/5)**

Call us : [800-304-5357](tel:800-304-5357)

## Product attributes

Antibody number	#2827
Antibody reactivity (target)	Cig2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	3A11/5
Isotype	IgG1, kappa
Molecular weight	~50 kDa
Synonyms	B type S phase cyclin antibody
Human gene symbol	Not Known
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Bacterial <i>S. pombe</i> Cig2 Purified protein.
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	<i>S. pombe</i>
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Red blood cell lysate; Human colon and lung tissue.
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	<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 11, 2025 at 11:36:40 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cig2-monoclonal-mouse-antibody-3a11-5/>