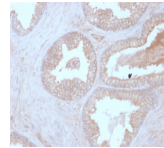


# Calcineurin B / PPP3R1 Monoclonal Mouse Antibody (CALNB/2342)



## Product Description

This antibody recognizes the Calcineurin B alpha protein, encoded by the PPP3R1 (Protein Phosphatase 3 Regulatory Subunit B Alpha) gene. Calcineurin is an enzyme that dephosphorylates serine and threonine residues in proteins. It is a heterodimer of a 59 kDa catalytic A subunit and a 19 kDa regulatory B subunit that is activated by the binding of calcium ions and calmodulin. Calcineurin is expressed in many tissues, but its levels are highest in the brain, where it may play a role in learning and memory. It has many substrates, including NFAT, a transcription factor that is activated by dephosphorylation. Complexes of the immuno-suppressants cyclosporine and FK506 with immunophilin proteins such as cyclophilin and FKBP12 are potent and specific inhibitors of Calcineurin activity. Alterations in Calcineurin activity are suspected to play a role in cardiac hypertrophy and graft versus host disease in organ transplantation. 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 2342, Calcineurin B / PPP3R1 Monoclonal Mouse Antibody (CALNB/2342)**

## Product attributes

Antibody number	#2342
Antibody reactivity (target)	Calcineurin B, PPP3R1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CALNB/2342
Isotype	IgG1, kappa
Molecular weight	19 kDa
Synonyms	Calcineurin B; Calcineurin subunit B type 1; Protein Phosphatase 3 Regulatory Subunit B Alpha; CALNB1; Cna2; CNB; CNB1; PPP3R1
Human gene symbol	PPP3R1
Entrez gene ID	5534
SwissProt	P63098
Unigene	280604
Immunogen	Recombinant full-length Calcineurin B protein
Antibody target cellular localization	Nucleus & cytoplasm
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Positive control	HeLa, MCF-7 or A431 cells. Prostate, Thyroid or Pancreas.
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 research areas	Ion channels
Cell/tissue expression	Brain, T-cells

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>
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.