Creatine Kinase B Monoclonal Mouse Antibody (CPTC-CKB-2)



Product Description

Creatine kinases (CK) are a large family of isoenzymes that regulate levels of ATP in subcellular compartments, where they provide ATP at sites of fluctuating energy demand by the transfer of phosphates between creatine and adenine nucleotides. CKs provide the energy of phosphate hydrolysis necessary to drive the normal function of many cellular systems. In cells, the cytosolic CK enzymes consist of two subunits, which can be either B (brain type) or M (muscle type). There are three different isoenzymes: CKMM, CKBB and CKMB. This MAb recognizes the CKBB isoenzyme and does not react with the B subunit in CKMB.

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 2233, Anti-Creatine Phosphokinase-BB (CPTC-CKB-2)

Product attributes

Product attributes				
Antibody number	#2233			
Antibody reactivity (target)	Creatine Phosphokinase-BB			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	CPTC-CKB-2			
Isotype	IgG2b, kappa			
Molecular weight	42 kDa			
Synonyms	BCK; Brain creatine kinase; Ckb; Creatine kinase B chain; Creatine kinase B-type; Creatine Kinase BB Isoenzyme; Creatine phosphokinase BB; Epididymis luminal protein 211; Epididymis secretory protein Li 29; HEL 211			
Human gene symbol	CKB			
Entrez gene ID	1152			
SwissProt	P12277			
Unigene	173724			
Immunogen	Recombinant human full-length CKB protein			
Immunogen Verified antibody applications	Recombinant human full-length CKB protein WB (verified)			
Verified antibody	y ,			
Verified antibody applications Antibody target cellular	WB (verified)			
Verified antibody applications Antibody target cellular localization	WB (verified) Cytoplasmic			
Verified antibody applications Antibody target cellular localization Species reactivity	WB (verified) Cytoplasmic Human			
Verified antibody applications Antibody target cellular localization Species reactivity Positive control	WB (verified) Cytoplasmic Human SH-SY5Y, HEK-293 and Cerebellum.			
Verified antibody applications Antibody target cellular localization Species reactivity Positive control Shipping condition	WB (verified) Cytoplasmic Human SH-SY5Y, HEK-293 and Cerebellum. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Verified antibody applications Antibody target cellular localization Species reactivity Positive control Shipping condition Storage Conditions	WB (verified) Cytoplasmic Human SH-SY5Y, HEK-293 and Cerebellum. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C			
Verified antibody applications Antibody target cellular localization Species reactivity Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	WB (verified) Cytoplasmic Human SH-SY5Y, HEK-293 and Cerebellum. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: BSA-free: 1 mg/mL in PBS/0.05% azide, Purified: BSA-free: 1 mg/mL in			

Chinese hamster ovary cells. Inquire for the specific lot.

Call us: 800-304-5357 Email: btinfo@biotium.com

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 November 3, 2025 at 11:09:18 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/creatine-kinase-b-monoclonal-mouse-antibody-cptc-ckb-2/