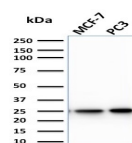


Calcineurin B homologous protein 2 / HCC Antigen 520 Monoclonal Mouse Antibody (CPTC-CHP2-1)



Product Description

Calcineurin B homologous protein 2, also known as hepatocellular carcinoma-associated antigen 520, is a 196 amino acid protein that contains 4 EF-hand domains and plays a potential role in transmembrane Na/H exchange. By binding to and activating NHE-1, CHP2 increases the pH and works to protect cells from serum deprivation-induced death. Though not typically detected in normal tissues, CHP2 is highly expressed in malignantly transformed cells and is therefore considered to be a tumor-associated antigen. Ectopic expression of CHP2 promotes proliferation of HEK293 cells and knockdown of CHP2 mRNA in HepG2 cells inhibits cell proliferation. Like calcineurin B, CHP2 can bind to and stimulate phosphatase activity of calcineurin A and activate the calcineurin/NFAT signaling pathway.

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 2266, Calcineurin B homologous protein 2 / HCC Antigen 520 Monoclonal Mouse Antibody (CPTC-CHP2-1)

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

Product attributes

Antibody number	#2266
Antibody reactivity (target)	Calcineurin B homologous protein 2, HCC Antigen 520
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-CHP2-1
Isotype	IgG2c, kappa
Molecular weight	22 kDa
Synonyms	Calcineurin B homologous protein 2; Calcineurin-like EF hand protein 2; HCA520; hepatocellular carcinoma antigen gene 520
Human gene symbol	CHP2
Entrez gene ID	63928
SwissProt	O43745
Unigene	178589
Immunogen	Recombinant human full-length CHP2 protein
Antibody target cellular localization	Nucleus & cytoplasm
Verified antibody applications	IF (verified), WB (verified)
Species reactivity	Human
Antibody application notes	Western blot 1-2 ug/mL; Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined.
Positive control	MCF-7, PC-3, HT29, K562 or Jurkat cell lysates. Human appendix or testis.
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
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	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 December 6, 2025 at 03:29:42 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/calcineurin-b-homologous-protein-2-hcc-antigen-520-monoclonal-mouse-antibody-cptc-chp2-1/>