Carbonic Anhydrase VIII Monoclonal Mouse Antibody (CPTC-CA8-2)



Product Description

Carbonic anhydrase VIII (carbonic anhydrase 8) was named because of sequence similarity to other known carbonic anhydrase genes. However, the gene product lacks carbonic anhydrase activity (i. e. , the reversible hydration of carbon dioxide). The gene product continues to carry a carbonic anhydrase designation based on clear sequence identity to other members of the carbonic anhydrase gene family. Mutations in the gene cause neurological disorders in mice.

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 2232, Carbonic Anhydrase VIII Monoclonal Mouse Antibody (CPTC-CA8-2)

Product attributes

Product attributes				
Antibody number	#2232			
Antibody reactivity (target)	•			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	CPTC-CA8-2			
Isotype	IgG2a, kappa			
Molecular weight	32.97 kDa			
Synonyms	Carbonic Anhydrase VIII; Carbonic Anhydrase 8; CA-VIII; CA-Related Protein			
Human gene symbol	CA8			
Entrez gene ID	767			
SwissProt	P35219			
Unigene	654388			
Immunogen	Recombinant full-length human CA8 protein			
Verified antibody applications	IHC (FFPE) (verified), WB (verified)			
Antibody target cellular localization	Cytoplasmic			
Species reactivity	Human			
Positive control	Human cerebellum. Selective expression in purkinje cells. HEK-293 and K-562 cell lines.			
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 (Bos taurus), or recombinant BSA produced 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 01:20:45 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/carbonic-anhydrase-viii-monoclonal-mouse-antibody-cptc-ca8-2/