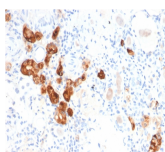


Calbindin 1 Monoclonal Mouse Antibody (CALB1/2782)



Product Description

Calbindin 1, also known as calbindin D28K, is a member of a family of calcium binding proteins. This protein family also includes calmodulin, calbindin 2, S-100, and troponin C. Calbindin 1 is a 261-amino acid protein with 6 EF-hand domains, 4 of which are active calcium-binding domains. Expressed in brain, ovary, uterus, testis, pancreas, liver, kidney and intestine, calbindin 1 acts as a calcium-buffering agent and alters the activity of the plasma membrane ATPase. In neuronal cells, calbindin 1 modulates calcium channel activity, calcium transients and intrinsic neuronal firing activity. Also, calbindin 1 has been implicated to play a role in apoptosis and microtubule function.

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 2782, Calbindin 1 Monoclonal Mouse Antibody (CALB1/2782)

Product attributes

Antibody number	#2782
Antibody reactivity (target)	Calbindin 1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CALB1/2782
Isotype	IgG2b, kappa
Molecular weight	52 kDa
Synonyms	Calbindin; Calbindin 1; Calbindin D28; Calbindin D28K; CALB1; Vitamin D dependent calcium binding protein
Human gene symbol	CALB1
Entrez gene ID	793
SwissProt	P05937
Unigene	65425
Immunogen	Recombinant fragment (around aa 7-96) of human CALB1 protein (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Human
Positive control	293T or HepG2 cells. Prostate, Lung, Pancreas, Kidney and Liver.
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
Validated in protein array	Monospecific
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 (<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	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 25, 2024 at 08:14:03 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/calbindin-1-monoclonal-mouse-antibody-calb1-2782/>