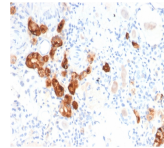


Calbindin 1 Monoclonal Mouse Antibody (CALB1/2782)



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 assay	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Antibody research areas	Ion channels

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
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.