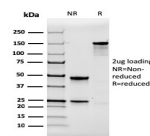


Calbindin 1 Monoclonal Mouse Antibody (CALB1/2364)



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 2364, Calbindin 1 Monoclonal Mouse Antibody (CALB1/2364)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #2364 |
| Antibody reactivity (target) | Calbindin 1 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | CALB1/2364 |
| Isotype | IgG1, 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) |
| Antibody target cellular localization | Nucleus & cytoplasm |
| Species reactivity | Human |
| Antibody application notes | For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry |
| 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 (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 August 24, 2025 at 02:27:42 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/calbindin-1-monoclonal-mouse-antibody-calb1-2364/>