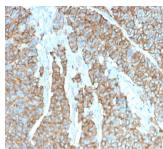


# Drebrin 1 Monoclonal Mouse Antibody (DBN1/3393)



## Product Description

Drebrins (developmentally regulated brain proteins) are cytoplasmic proteins that bind F-actin in the brain and are involved in cell migration, extension of neuronal processes and plasticity of dendrites. There are three isoforms: two embryonic types (E1 and E2); and an adult type (A), generated by alternative RNA splicing from a single Drebrin gene. Drebrins are expressed mainly in brain neurons but are also found in skeletal muscle, heart, placenta, pancreas and kidney. Drebrin has been designated as a marker of the dendritic spine. Decreases in Drebrin levels in the brain have been associated with Alzheimer's disease and Down syndrome.

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 3393, Anti-Drebrin 1 (DBN1/3393)**

## Product attributes

|                                       |   |
|---------------------------------------|---|
| Antibody number                       | #3393   |
| Antibody reactivity (target)          | Drebrin 1   |
| Antibody type                         | Primary   |
| Host species                          | Mouse   |
| Clonality                             | Monoclonal  |
| Clone                                 | DBN1/3393   |
| Isotype                               | IgG1, kappa   |
| Molecular weight                      | 120 kDa   |
| Synonyms                              | Developmentally regulated brain protein; Drebrin 1; Drebrin; Drebrin E; Drebrin E2; Drebrin1; DrebrinE  |
| Human gene symbol                     | DBN1  |
| Entrez gene ID                        | 1627  |
| SwissProt                             | Q16643  |
| Unigene                               | 130316  |
| Immunogen                             | Recombinant fragment (aa 150-281) of human DBN1 protein (exact sequence is proprietary)   |
| Verified antibody applications        | IHC (FFPE) (verified), WB (verified)  |
| Antibody target cellular localization | Cytoskeleton, Plasma membrane   |
| Species reactivity                    | Human   |
| Positive control                      | Brain neurons. Also found in the heart, placenta, skeletal muscle, kidney and pancreas.   |
| 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   |
| 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.                   |
| Cell/tissue expression                | Heart, Skeletal Muscle, Brain, Kidney, Neurons, Pancreas, Placenta  |

| Antibody # prefix | Conjugation        | Ex/Em (nm) | Laser line | Detection channel        | Dye Features                     |
|-------------------|--------------------|------------|------------|--------------------------|----------------------------------|
| BNC04             | CF®405S            | 404/431    | 405        | DAPI (microscopy), AF405 | <a href="#">CF®405S Features</a> |
| BNC88             | CF®488A            | 490/515    | 488        | GFP, FITC                | <a href="#">CF®488A Features</a> |
| BNC68             | CF®568             | 562/583    | 532, 561   | RFP, TRITC               | <a href="#">CF®568 Features</a>  |
| BNC94             | CF®594             | 593/614    | 561        | Texas Red®               | <a href="#">CF®594 Features</a>  |
| BNC40             | CF®640R            | 642/662    | 633-640    | Cy®5                     | <a href="#">CF®640R Features</a> |
| BNC47             | CF®647             | 650/665    | 633-640    | Cy®5                     | <a href="#">CF®647 Features</a>  |
| BNC74             | CF®740             | 742/767    | 633-685    | 775/50                   | <a href="#">CF®740 Features</a>  |
| 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 July 18, 2025 at 05:07:38 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/drebrin-1-monoclonal-mouse-antibody-dbn1-3393/>