## FABP5 Monoclonal Mouse Antibody (CPTC-FABP5-3)



## **Product Description**

This gene encodes the fatty acid binding protein found in epidermal cells, and was first identified as being upregulated in psoriasis tissue. Fatty acid binding proteins are a family of small, highly conserved, cytoplasmic proteins that bind long-chain fatty acids and other hydrophobic ligands. FABPs may play roles in fatty acid uptake, transport, and metabolism. Polymorphisms in this gene are associated with type 2 diabetes. The human genome contains many pseudogenes similar to this locus. FABP5 is also associated with poor survival in triple-negative breast cancer. Additionally, FABP5 gene is upregulated in colorectal cancer cells compared to normal colon cells in a manner that correlates with disease stage and that FABP5 significantly promotes colorectal cancer cell growth and metastatic potential.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2225, Anti-FABP5 (CPTC-FABP5-3)

## Product attributes

| Product attributes                       |   |  |  |
|--|---|--|--|
| Antibody number                          | #2225   |  |  |
| Antibody reactivity (target)             | FABP5   |  |  |
| Antibody type                            | Primary   |  |  |
| Host species                             | Mouse   |  |  |
| Clonality                                | Monoclonal  |  |  |
| Clone                                    | CPTC-FABP5-3  |  |  |
| Isotype                                  | IgG2a, kappa  |  |  |
| Molecular weight                         | 15.4 kDa  |  |  |
| Synonyms                                 | FABP5; E-FABP; EFABP; PA-FABP; PAFABP   |  |  |
| Human gene symbol                        | FABP5   |  |  |
| Entrez gene ID                           | 2171  |  |  |
| SwissProt                                | Q01469  |  |  |
| Unigene                                  | 408061  |  |  |
| Immunogen                                | Recombinant human full-length FABP5 protein   |  |  |
| Verified antibody applications           | IHC (FFPE) (verified), WB (verified)  |  |  |
| Antibody target cellular<br>localization | Nucleus & cytoplasm   |  |  |
| Species reactivity                       | Human   |  |  |
| Positive control                         | MOLT-4 and HEK-293 cell lines. Esophagus, Skin.<br>Keratinocytes; highly expressed in psoriatic skin. Colon.  |  |  |
| Shipping condition                       | Room temperature  |  |  |
| Storage Conditions                       | Store at 2 to 8 $^{\circ}$ C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ 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),<br>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,<br>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 4, 2025 at 08:04:32 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/fabp5-monoclonal-mouse-antibody-cptc-fabp5-3/">https://biotium.com/product/fabp5-monoclonal-mouse-antibody-cptc-fabp5-3/</a>