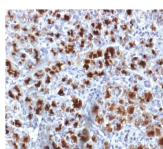


# CELA3B / ELA3B Monoclonal Mouse Antibody (CELA3B/1757)



## Product Description

This MAbs recognizes a protein of ~30 kDa, identified as CELA3B (Chymotrypsin like elastase family member 3B). Elastases form a subfamily of serine proteases that hydrolyze many proteins in addition to elastin. Humans have six elastase genes which encode the structurally similar proteins elastase 1, 2, 2A, 2B, 3A, and 3B. Unlike other elastases, elastase 3B has little elastolytic activity. Like most of the human elastases, elastase 3B is secreted from the pancreas as a zymogen and, like other serine proteases such as trypsin, chymotrypsin and kallikrein; it has a digestive function in the intestine. Elastase 3B preferentially cleaves proteins after alanine residues. Elastase 3B may also function in the intestinal transport and metabolism of cholesterol. Both elastase 3A and elastase 3B have been referred to as protease E and as elastase 1, and excretion of this protein in fecal material is frequently used as a measure of pancreatic 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 1757, Anti-CELA3B (CELA3B/1757)**

## Product attributes

|  |   |
|--|---|
| <b>Antibody number</b>                       | #1757   |
| <b>Antibody reactivity (target)</b>          | CELA3B, ELA3B   |
| <b>Antibody type</b>                         | Primary   |
| <b>Host species</b>                          | Mouse   |
| <b>Clonality</b>                             | Monoclonal  |
| <b>Clone</b>                                 | CELA3B/1757   |
| <b>Isotype</b>                               | IgG1, kappa   |
| <b>Molecular weight of antigen</b>           | 17 kDa  |
| <b>Synonyms</b>                              | Chymotrypsin like elastase family member 3B (CELA3B); ELA3B; Elastase IIIb; Protease E  |
| <b>Human gene symbol</b>                     | CELA3B  |
| <b>Entrez gene ID</b>                        | 23436   |
| <b>SwissProt</b>                             | P08861  |
| <b>Unigene</b>                               | 728752  |
| <b>Immunogen</b>                             | Recombinant human CELA3B protein fragment (aa82-238) (exact sequence is proprietary)  |
| <b>Verified antibody applications</b>        | IHC (FFPE) (verified)   |
| <b>Antibody target cellular localization</b> | Cytoplasmic   |
| <b>Species reactivity</b>                    | Human   |
| <b>Antibody application notes</b>            | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, ELISA: 2-4 ug/mL for coating order Ab without BSA, Immunohistology (formalin) 1-2 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 for 10-20 min followed by cooling at RT for 20 min, Immunofluorescence 0.5-1 ug/mL Western blotting 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user |
| <b>Positive control</b>                      | Panc-28 or PANC1 cells. Pancreas  |
| <b>Shipping condition</b>                    | Room temperature  |
| <b>Storage Conditions</b>                    | Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C   |
| <b>Shelf life</b>                            | Guaranteed for at least 24 months from date of receipt when stored as recommended   |
| <b>Regulatory status</b>                     | For research use only (RUO)   |
| <b>Antibody/conjugate formulation</b>        | 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   |
| <b>Validated in protein array</b>            | Monospecific  |
| <b>Antibody research areas</b>               | Proteases   |
| <b>Product origin</b>                        | 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 | <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 June 10, 2026 at 01:58:19 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/mouse-monoclonal-anti-cela3b-ela3b-cela3b1757/>