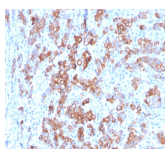


# TIM3 / HAVCR2 / CD366 Monoclonal Mouse Antibody (TIM3/3113)



## Product Description

TIMs are type I transmembrane glycoproteins with one Ig-like V-type domain and a Ser/Thr-rich mucin stalk. TIM-3 is expressed on the surface of effector T cells (CD4 Th1 and CD8 Tc1) but not on helper T cells (CD4 Th2 and CD8 Tc2). In chronic inflammation, autoimmune disorders, and some cancers, TIM-3 is upregulated on several other hematopoietic cell types. The Ig domain of TIM-3 interacts with a ligand on resting but not activated Th1 and Th2 cells. The glycosylated Ig domain of TIM-3 binds cell-associated galectin-9. This induces TIM-3 Tyr phosphorylation and pro-apoptotic signaling. TIM-3 functions as a negative regulator of Th1 cell activity. Its blockade results in increased IFN-gamma production, Th1 cell proliferation and cytotoxicity, regulatory T cell development, and increases in macrophage and neutrophil infiltration into sites of inflammation.

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 3113, Anti-CD366|HAVCR2|TIM3 (TIM3/3113)**

## Product attributes

|                                       |  |
|---------------------------------------|--|
| Antibody number                       | #3113  |
| Antibody reactivity (target)          | CD366, HAVCR2, TIM3  |
| Antibody type                         | Primary  |
| Host species                          | Mouse  |
| Clonality                             | Monoclonal   |
| Clone                                 | TIM3/3113  |
| Isotype                               | IgG2a, kappa   |
| Molecular weight                      | 60 kDa   |
| Synonyms                              | CD366; HAVR2; Hepatitis A virus cellular receptor 2 (HAVCR2); Kidney injury molecule 3 (KIM3); T-cell immunoglobulin and mucin domain-containing protein 3; T-cell immunoglobulin mucin receptor 3; T-cell membrane protein 3; TIM3; TIMD3 |
| Human gene symbol                     | HAVCR2   |
| Entrez gene ID                        | 84868  |
| SwissProt                             | Q8TDQ0   |
| Unigene                               | 710500   |
| Immunogen                             | Recombinant fragment of human TIM3 protein (around aa 22-202) (exact sequence is proprietary)  |
| Verified antibody applications        | IHC (FFPE) (verified)  |
| Antibody target cellular localization | Plasma membrane  |
| Species reactivity                    | Human  |
| Positive control                      | PC3, BT474, HepG2, HDLM-2 or Daudi cells. Tonsil, lymph node or spleen.  |
| 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.  |

| 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 August 7, 2025 at 07:32:40 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/tim3-havcr2-cd366-monoclonal-mouse-antibody-tim3-3113/>