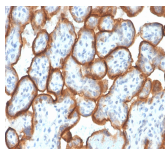


## VISTA / Gi24 Monoclonal Mouse Antibody (VISTA/3007)



### Product Description

VISTA / Gi24 is a transmembrane protein expressed in bone, on embryonic stem cells (ESCs), and on tumor cell surfaces. On ESC s, Gi24 appears to positively interact with BMP-4, potentiating BMP signaling and the transition from an undifferentiated to a differentiated state. On tumor cells, Gi24 both promotes MT1-MMP expression and activity and serves as a substrate for MT1-MMP. This increases the potential for cell motility. Mature human Gi24 contains a 162aa extracellular region with one V-type Ig-like domain and a 96aa cytoplasmic domain. Human Gi24 undergoes proteolytic cleavage by MT1-MMP, generating a soluble 30 kDa extracellular fragment plus a 25-30 kDa membrane-bound fragment. VISTA is a negative checkpoint regulator and is expressed on myeloid cells, T-cells and human TiLs (tumor infiltrating lymphocytes) on MDSCs (myeloid-derived suppressor cells) in the TME (tumor microenvironment). It is very likely both a ligand and receptor and is a promising target for cancer immunotherapy.

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 3007, Anti-B7-H5|VISTA|VSIR (VISTA/3007)**

### Product attributes

|                                       |  |
|---------------------------------------|--|
| Antibody number                       | #3007  |
| Antibody reactivity (target)          | B7-H5, VISTA, VSIR   |
| Antibody type                         | Primary  |
| Host species                          | Mouse  |
| Clonality                             | Monoclonal   |
| Clone                                 | VISTA/3007   |
| Isotype                               | IgG2b, kappa   |
| Molecular weight                      | 34 kDa   |
| Synonyms                              | B7H5; DD1alpha; Gi24; PD-1H; PDCD1 homolog; Platelet receptor Gi24; PP2135; Stress-induced secreted protein 1 (SISP1); V domain Ig suppressor of T cell activation (VISTA); V Set Immuno-Regulatory Receptor |
| Human gene symbol                     | VSIR   |
| Entrez gene ID                        | 64115  |
| SwissProt                             | Q9H7M9   |
| Unigene                               | 47382  |
| Immunogen                             | Recombinant full-length human VISTA protein  |
| Verified antibody applications        | IHC (FFPE) (verified)  |
| Antibody target cellular localization | Plasma membrane  |
| Species reactivity                    | Human  |
| Positive control                      | Spleen or Placenta.  |
| 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 December 13, 2025 at 10:55:56 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/vista-gi24-monoclonal-mouse-antibody-vista-3007/>