Glowing products for science

CD57 / B3GAT1 Monoclonal Mouse Antibody (NK-1)



Product Description

Anti-CD57 marks a subset of lymphocytes known as natural killer (NK) cells. Follicular center cell lymphomas often contain many NK cells within the neoplastic follicles. Anti-CD57 also stains neuroendocrine cells and their derived tumors, including carcinoid tumor and medulloblastoma. Anti-CD57 can also be useful in separating type B3 thymoma from thymic carcinoma when combined with a panel that includes antibodies against GLUT1, CD5, and CEA.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL. Catalog number key for antibody number 0016, Anti-B3GAT1|CD57 (NK-1)

Call us: 800-304-5357 Email: btinfo@biotium.com

Product attributes

| Antibody number | #0016 | | | |
|--|--|--|--|--|
| Antibody reactivity (target) | B3GAT1, CD57 | | | |
| Antibody type | Primary | | | |
| Host species | Mouse | | | |
| Clonality | Monoclonal | | | |
| Clone | NK-1 | | | |
| Isotype | IgM, kappa | | | |
| Molecular weight | ~110 kDa (Glycoprotein) | | | |
| Synonyms | 3-Glucuronyltransferase 1; B3GAT1; Galactosylgalactosylxylosylprotein 3-beta-Glucuronosyltransferase 1; GLCATP; GlcUAT-P; Glucuronosyltransferase P; UDP GlcUA Glycoprotein beta 1, 3 Glucuronyltransferase | | | |
| Human gene symbol | B3GAT1 | | | |
| Entrez gene ID | 27087 | | | |
| SwissProt | Q9P2W7 | | | |
| Unigene | 381050 | | | |
| Immunogen | Human peripheral blood mononuclear cells | | | |
| Verified antibody applications | B3GAT1 27087 Q9P2W7 381050 Human peripheral blood mononuclear cells IHC (FFPE) (verified) Plasma membrane Human | | | |
| Antibody target cellular localization | Plasma membrane | | | |
| Species reactivity | Human | | | |
| Expected antibody applications | Flow, surface (published for clone), IF (published for clone), WB (published for clone) | | | |
| Antibody application notes | primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Does not react with rat; others not known, Immunohistology formalin-fixed 2-4 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, | | | |
| Positive control | Lymph node or Tonsil | | | |
| Shipping condition | Room temperature | | | |
| Storage Conditions | Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C | | | |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended | | | |
| 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 | | | |
| Antibody research areas | Cancer, Immunology | | | |
| 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 | 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, 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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

- Asian J Androl (2010) 12(4): 548-555. (WB blot; Flow)
 Cancer Immunol Immunother (2011) 60:1683-1695. (Flow, surface)
 Cancer Sci (2016) 107: 846-852. (IHC, FFPE)
 Cell Stem Cell (2017) 20(6): 874-890.e7. (IF; Flow, surface)

This datasheet was generated on November 8, 2025 at 07:23:52 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-cd57-b3gat1-nk-1/