

CD16 / Fc Gamma Receptor III Monoclonal Mouse Antibody (HuNK2)

Product Description

Recognizes a protein of 50-65 kDa, identified as CD16 (Workshop IV; Code N39) (also known low affinity Fc receptor III for IgG (FcRIII) or Leu 11). CD16 exists as a polypeptide-anchored form (FCRIIIA or CD16A) on human natural killer (NK) cells and monocytes/ macrophages and as a glycosylphosphatidylinositol (GPI)-anchored form (FcRIIIB or CD16B) on neutrophils. CD16B is polymorphic and the two alleles are termed NA1 and NA2.3 CD16 plays a role in signal transduction, NK cell activation and antibody-dependent cellular cytotoxicity. This MAb has been showed to inhibit the binding of immune complex to NK cells, inhibit cytotoxicity of NK cells, and induce calcium fluxes in NK cells and neutrophils.

Also view our [CD16 Mouse Monoclonal Biotium Choice Antibody \(3G8\)](#). The antibody belongs to the Biotium Choice list of select antibodies that have been developed and validated in-house for optimal performance. The antibody has been validated in flow cytometry.

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 to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0928, Anti-CD16|Fc gamma receptor III (HuNK2)

Product attributes

| | |
|--|--|
| Antibody number | #0928 |
| Antibody reactivity (target) | CD16, Fc gamma receptor III |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | HuNK2 |
| Isotype | IgG2a, kappa |
| Molecular weight of antigen | 50-80 kDa |
| Synonyms | CD162; CLA; Cutaneous lymphocyte associated antigen; P-selectin glycoprotein ligand 1; PSGL1; Selectin P ligand; SELPLG |
| Human gene symbol | FCGR3A |
| Entrez gene ID | 2214 |
| SwissProt | P08637 |
| Unigene | 372679 |
| Immunogen | Human peripheral blood lymphocytes from a NK-leukemia patient. |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Expected antibody applications | Flow, surface (published for clone) |
| Antibody application notes | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 0.5-1 ug/mL, Functional Studies order Ab without BSA & azide). Optimal dilution for a specific application should be determined by user |
| Positive control | K-562 or U-937 cells. Lymph nodes and tonsils. |
| 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 |
| 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 | 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. |
| Cell/tissue expression | Monocytes/macrophages, Neutrophils, NK cells |

| 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.

- Blood (2001) 98(1):140-145. (Flow, surface)
- Blood (2002) 100(13): 4512-4520. (cell depletion)