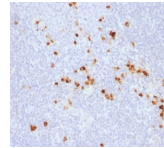


Human IgG Heavy Chain Recombinant Monoclonal Rabbit Antibody (IG507R)



Product Description

This antibody recognizes a protein of 75 kDa, identified as gamma heavy chain of human immunoglobulins. It does not cross-react with alpha (IgA), mu (IgM), epsilon (IgE), or delta (IgD), heavy chains, T-cells, monocytes, granulocytes, or erythrocytes. This antibody is useful in the identification of leukemias, plasmacytomas, and certain non-Hodgkin's lymphomas. The most common feature of these malignancies is the restricted expression of a single heavy chain class. Demonstration of clonality in lymphoid infiltrates indicates that the infiltrate is clonal and therefore malignant. 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.

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #0507 |
| Antibody reactivity (target) | Human IgG Heavy Chain |
| Antibody type | Primary, Anti-Human Immunoglobulin |
| Host species | Rabbit |
| Clonality | Recombinant Monoclonal |
| Clone | IG507R |
| Isotype | IgG, kappa |
| Molecular weight | 75 kDa |
| Synonyms | G1m Marker; G2m Marker; G3m Marker; G4m Marker; HDC; Heavy Chain Disease Protein; Human Immunoglobulin G; Ig gamma1/2/3/4 Chain C Region;IGHG1; IGHG2; IGHG3; IGHG4; Immunoglobulin Heavy Constant 1/2/3/4 |
| Human gene symbol | IGHG |
| Entrez gene ID | 3500 (IGHG1), 3501 (IGHG2), 3502 (IGHG3), 3503 (IGHG4) (Human) |
| SwissProt | P01857, P01859, P01860, P01861 |
| Unigene | 510635 |
| Immunogen | Purified human Ig Gamma Chain |
| Antibody target cellular localization | Secreted (extracellular) |
| Verified antibody applications | IHC (FFPE) (verified) |
| Species reactivity | Human |
| 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, Immunohistochemistry (formalin): 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user |
| Positive control | 293T, Raji or hPBL cells. Tonsil 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 |
| 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 |
| Validated in protein | Monospecific |
| Antibody research areas | Cancer, Immunology, Inflammation |
| Cell/tissue expression | B-cells |
| Tumor expression | Leukemia/lymphoma |

| Antibody # prefix | Conjugation | Ex/Em (nm) | Laser line | Detection channel |
|-------------------|--------------------|------------|------------|--------------------------|
| BNC04 | CF®405S | 404/431 | 405 | DAPI (microscopy), AF405 |
| BNC88 | CF®488A | 490/515 | 488 | GFP, FITC |
| BNC68 | CF®568 | 562/583 | 532, 561 | RFP, TRITC |
| BNC94 | CF®594 | 593/614 | 561 | Texas Red® |
| BNC40 | CF®640R | 642/662 | 633-640 | Cy®5 |
| BNC47 | CF®647 | 650/665 | 633-640 | Cy®5 |
| 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 |

Dye Features

- [CF®405S Features](#)
- [CF®488A Features](#)
- [CF®568 Features](#)
- [CF®594 Features](#)
- [CF®640R Features](#)
- [CF®647 Features](#)

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, LI-COR Bioscience.