



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: [btinfo@biotium.com](mailto:btinfo@biotium.com)

## Human IgG Heavy Chain Monoclonal Mouse Antibody (IG217 + IG266)

### Product Description

This MAb is specific to heavy chain of IgG and shows minimal cross-reaction with heavy chains of other immunoglobulins. It is reactive with all subclasses of Gamma heavy chain(IgG1, IgG2a, IgG2b, IgG3). The most common feature of plasmacytomas, and certain non-Hodgkin's lymphomas 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. Ab-3 is superb for staining of formalin-fixed tissues.

#### Product attributes

Antibody number	#1262
Antibody type	Primary, Anti-Human Immunoglobulin
Host species	Mouse
Clonality	Monoclonal
Clone	IG217 + IG266
Isotype	IgG1, kappa + IgG2a, 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; 3501; 3502; 3503
SwissProt	P01857, P01859, P01860, P01861
Unigene	510635
Immunogen	Purified human IgG
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Antibody application notes	Flow cytometry: 1 ug/million cells. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunohistochemistry (formalin-fixed): 0.5-1 ug/mL for 30 minutes at RT, Immunofluorescence: 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 minutes followed by cooling at RT for 20 minutes, 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
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Cell/tissue expression	B-cells
Tumor expression	Leukemia/lymphoma
Antibody research areas	Cancer, Immunology, Inflammation

This datasheet was generated on April 25, 2024 at 02:48:04 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/human-igg-heavy-chain-monoclonal-mouse-antibody-ig217-ig266/>