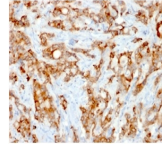


Beta-2 Microglobulin Recombinant Monoclonal Mouse Antibody (rB2M/961)



Product Description

Beta-2 microglobulin is a component of the class I major histocompatibility complex (MHC I) where it functions in antigen processing and presentation. Involved in the presentation of peptide antigens to the immune system. Human Beta-2 microglobulin is expressed on many types of cells including lymphocytes, thymocytes, monocytes, granulocytes, platelets, endothelial cells, and epithelial cells. It is absent on erythrocytes. This MAb is specific to human Beta-2 microglobulin and does not react with non-human primate cells. This antibody reacts with all cell types excluding erythrocytes. The detection of Beta-2 microglobulin in body fluids has been used as a tumor marker and for monitoring patients with HIV infection. 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 1537, Anti-Beta-2 Microglobulin (rB2M/961)**

Product attributes

Antibody number	#1537
Antibody reactivity (target)	Beta-2 Microglobulin
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rB2M/961
Isotype	IgG2b, kappa
Molecular weight	12 kDa
Synonyms	B2M; Beta 2 microglobulin
Human gene symbol	B2M
Entrez gene ID	567
SwissProt	P61769
Unigene	534255
Immunogen	Recombinant human full-length B2M protein
Antibody target cellular localization	Endoplasmic reticulum, Plasma membrane, Secreted (extracellular)
Verified antibody applications	Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified), WB (verified)
Species reactivity	Human, Non-human primates
Positive control	HL-60 or HeLa cells. Melanomas and Lymphoma. Carcinoma of Stomach, Cervix, Endometrial, Kidney or Colon.
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 array	Monospecific
Antibody research areas	Cancer, Immunology, Inflammation

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
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, LI-COR Bioscience.