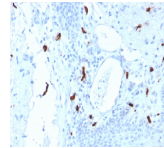


Tryptase Monoclonal Mouse Antibody (TPSAB1/1961)



Product Description

Tryptases comprise a family of trypsin-like serine proteases (peptidase family S1). Tryptases are stored in mast cell secretory granules and basophils. Mast cells are connective tissue cells derived from blood-forming tissues that line arterial walls and secrete substances, which mediate inflammatory and immune responses. Tryptases are released into the extracellular environment and are resistant to all known endogenous proteinase inhibitors. This antibody reacts with mast cells distributed in skin, synovium, lung, and heart. This antibody does not bind with any other cell type. Human mast cell tryptase is considered to be an important marker of mast cell activation and is an important mediator of inflammation. Mastocytosis is a term collectively used for a group of disorders in which there is abnormal accumulation of mast cells in one or multiple organs. Anti-tryptase, combined with anti-CD2, anti-CD25, and anti-CD117, can be useful in identifying reactive mast cell hyperplasia, myelogenous neoplasms, mast cell leukemia, and mastocytosis.

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 1961, Anti-Tryptase (TPSAB1/1961)

Product attributes

Antibody number	#1961
Antibody reactivity (target)	Tryptase
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	TPSAB1/1961
Isotype	IgG1, kappa
Molecular weight	31-36 kDa
Synonyms	Lung Tryptase; Mast cell alpha II Tryptase; Mast cell beta I Tryptase; Mast cell protease 7; Mast cell protease II; MCP7; Pituitary Tryptase; Skin Tryptase; TPS1; TPSAB1; TPSB1; Tryptase 1; Tryptase alpha 1; Tryptase skin
Human gene symbol	TPSAB1
Entrez gene ID	7177
SwissProt	Q15661
Unigene	405479
Immunogen	Recombinant human Tryptase protein fragment (around aa 115-233) (Exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Positive control	Skin or Tonsil
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	Immunology, Inflammation
Tumor expression	Leukemia/lymphoma

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, and LI-COR Bioscience.

This datasheet was generated on September 2, 2024 at 12:21:57 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/tryptase-monoclonal-mouse-antibody-tpsab1-1961/>