HA Tag Monoclonal Mouse Antibody (16.43)



Product Description

Human influenza hemagglutinin (HA) is a surface glycoprotein required for the infectivity of the human virus. The HA tag is derived from the HA molecule corresponding to amino acids 98-106 has been extensively used as a general epitope tag in expression vectors. Many recombinant proteins have been engineered to express the HA tag, which does not appear to interfere with the bioactivity or the biodistribution of the recombinant protein. This tag facilitates the detection, isolation, and purification of the proteins

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 2543, Anti-HA Tag (16.43)

Product attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#2543			
Antibody reactivity (target)	A Tag			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	16.43			
Isotype	IgG1, kappa			
Molecular weight	60 kDa			
Synonyms	HA epitope tag; HA1; HA2; hemagglutinin; Hemagglutinin HA1 chain; Hemagglutinin HA2 chain			
Human gene symbol	Not Known			
Entrez gene ID	Not Known			
SwissProt	Not Known			
Unigene	Not Known			
Immunogen	Semi-purified mitochondrial preparation			
Verified antibody applications	WB (verified)			
Species reactivity	Human			
Species reactivity Antibody application notes	Human Does not react with Mouse or Rat; ELISA; Use Ab at 2-4 ug/mL for coating; Order Ab without BSA; Optimal dilution for a specific application should be determined by user., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody			
	Does not react with Mouse or Rat; ELISA: Use Ab at 2-4 ug/mL for coating; Order Ab without BSA; Optimal dilution for a specific application should be determined by user., Higher concentration may be required for direct detection using primary antibody			
Antibody application notes	Does not react with Mouse or Rat; ELISA: Use Ab at 2-4 ug/mL for coating; Order Ab without BSA; Optimal dilution for a specific application should be determined by user., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody			
Antibody application notes Positive control	Does not react with Mouse or Rat; ELISA: Use Ab at 2-4 ug/mL for coating; Order Ab without BSA; Optimal dilution for a specific application should be determined by user., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody HeLa or HepG2 cells. Hepatic carcinoma.			
Antibody application notes Positive control Shipping condition	Does not react with Mouse or Rat; ELISA: Use Ab at 2-4 ug/mL for coating; Order Ab without BSA; Optimal dilution for a specific application should be determined by user., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody HeLa or HepG2 cells. Hepatic carcinoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Antibody application notes Positive control Shipping condition Storage Conditions	Does not react with Mouse or Rat; ELISA: Use Ab at 2-4 ug/mL for coating; Order Ab without BSA; Optimal dilution for a specific application should be determined by user., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody HeLa or HepG2 cells. Hepatic carcinoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C			
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	Does not react with Mouse or Rat; ELISA: Use Ab at 2-4 ug/mL for coating; Order Ab without BSA; Optimal dilution for a specific application should be determined by user., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody HeLa or HepG2 cells. Hepatic carcinoma. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA/, Upflied: 0.2 mg/mL in PBS/0.05% BSA/, Upflied: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflied: 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA/0.05% azide, Purflied: 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA/0.05% azide, Purflied: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflied: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purflied: 0.2 mg/mL in PBS/0.05% BSA/0.05% BSA/			

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.

This datasheet was generated on November 29, 2025 at 08:25:14 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/ha-tag-monoclonal-mouse-antibody-16-43/