Alpha-Fetoprotein Monoclonal Mouse Antibody (C2 + C3 + MBS-12)



Product Description

Alpha-fetoprotein (AFP) is an oncofetal glycoprotein with a single chain of 70 kDa. This MAb is highly specific to AFP and shows no cross-reaction with other oncofetal antigens or serum albumin. AFP is normally synthesized in the liver, intestinal tract, and yolk sac of the fetus. Antibody to AFP has been shown to be useful in detecting hepatocellular carcinomas (HCC) and germ cell neoplasms, especially yolk sac tumors.

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 0810, Anti-AFP (C2 C3 MBS-12)

Call us: 800-304-5357

Product attributes			
research-areas	Cancer, Developmental biology		
Antibody number	#0810		
Antibody reactivity (target)	AFP, Alpha-fetoprotein		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	C2 C3 MBS-12		
Isotype	IgG's		
Molecular weight	70 kDa		
Synonyms	Alpha-fetoglobulin; FETA; HPAFP		
Human gene symbol	AFP		
Entrez gene ID	174		
SwissProt	P02771		
Unigene	518808		
Immunogen	Recombinant full-length human Alpha fetoprotein		
Antibody target cellular localization	Cytoplasmic, Golgi apparatus		
Verified antibody applications	Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified)		
Species reactivity	Human		
Antibody application notes	Immunohistology formalin-fixed 1-2 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 minutes, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/mllion cells/0.1 mL, Optimal dilution for a specific application should be determined by user		
Positive control	HepG2 cells. Fetal liver or hepatocellular carcinoma		
Shipping condition	Room temperature		
	Room temperature		
Storage Conditions	Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C		
	Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when		
Storage Conditions Shelf life	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended		
Storage Conditions Shelf life Regulatory status Antibody/conjugate	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purifled: 0.2 mg/mL in PBS/0.05% BSA, Purifled; 0.2 mg/mL in PBS/0.05% BSA, Purifled, BSA-free: 1 mg/mL in		
Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA-free: 1 mg/mL in PBS without azide		

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 30, 2025 at 02:54:40 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-afp-alpha-fetoprotein-afp810/