Fibrinogen Alpha Chain Monoclonal Mouse Antibody (UC45)



Product Description

The plasma glycoprotein fibrinogen is synthesized in the liver and comprises three structurally different subunits: . Fibrinogen is important in platelet aggregation, the final step of the coagulation cascade (i. e. the formation of fibrin) and determination of plasma viscosity and erythrocyte aggregation. It is both constitutively expressed and inducible during an acute phase reaction. Hemostasis following tissue injury deploys essential plasma procoagulants (prothrombin and factors X, IX, V and VIII), which are involved in a blood coagulation cascade leading to the formation of insoluble fibrin clots and the promotion of platelet aggregation. Following vascular injury, fibrinogen is cleaved by thrombin to form fibrin, which is the most abundant component of blood clots. The cleavage products of fibrinogen regulate cell adhesion and spreading, display vasoconstrictor and chemotactic activities, and are mitogens for several cell types.

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 2849, Anti-Fibrinogen Alpha Chain (UC45)

Product attributes

Call us: 800-304-5357

Product attributes	
Antibody number	#2849
Antibody reactivity (target)	Fibrinogen Alpha Chain
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	UC45
Isotype	IgM, kappa
Molecular weight	~30 kDa
Synonyms	Ac1873; Fba5e; FGA; Fib; Fib2; Fibrinogen alpha chain; Fibrinogen; A alpha polypeptide; Fibrinogen-alpha polypeptide chain; Fibrinopeptide A
Human gene symbol	FGA
Entrez gene ID	2243
SwissProt	P02671
Unigene	351593
Immunogen	Human acute monoblastic leukemia cells.
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Expected antibody applications	Flow (published for clone), IHC (FFPE) (published for clone), IF (published for clone), WB (published for clone)
Antibody application notes	ELISA: For coating, order antibody without BSA; Flow cytometry: 0.5-1 ug/million cells; Immunofluorescence: 0.5-1 ug/mil.; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	Plasma.
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35°C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
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
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Email: btinfo@biotium.com

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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

- 1. J Exp Med (1983) 157: 473-485. (WB; IF)
- 2. Cancer Sci (2019) 110(8): 2658-2666. (IHC, FFPE)
- 3. Neurosci Lett (2015) 602:126-132. (IF; WB)
- 4. Artificial Cells, Nanomedicine, Biotechnology (2020) 48(1): 28-36. (Flow)

This datasheet was generated on November 4, 2025 at 10:49:10 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/fibrinogen-alpha-chain-monoclonal-mouse-antibody-uc45/