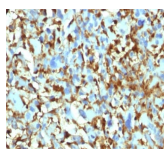


Factor XIIIa Monoclonal Mouse Antibody (F13A1/1448)



Product Description

This antibody recognizes a protein of 83 kDa, which is identified as Factor XIIIa. It has been identified in platelets, megakaryocytes, and fibroblast-like mesenchymal or histiocytic cells in the placenta, uterus, and prostate, monocytes and macrophages and dermal dendritic cells. Anti-factor XIIIa has been found to be useful in differentiating between dermatofibroma (almost all cases are positive), dermatofibrosarcoma protuberans (-/-) and desmoplastic malignant melanoma (-). Anti-factor XIIIa positivity is also seen in capillary hemangioblastoma, hemangioendothelioma, hemangiopericytoma, xanthogranuloma, xanthoma, hepatocellular carcinoma, glomus tumor, and meningioma.

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 1448, Anti-Factor-XIIIa (F13A1/1448)

Product attributes

Antibody number	#1448
Antibody reactivity (target)	Factor-XIIIa
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	F13A1/1448
Isotype	IgG2b, kappa
Molecular weight	83 kDa
Synonyms	Coagulation factor XIII A chain; Coagulation factor XIII A1 polypeptide; Coagulation factor XIIIa; F13A; F13a1; Factor XIIIa; Fibrin stabilizing factor, A subunit; Fibrinolytic; F5F, A subunit; Protein-glutamine gamma-glutamyltransferase A chain; TGase; Transglutaminase A chain; Transglutaminase. plasma
Human gene symbol	F13A1
Entrez gene ID	2162
SwissProt	P00488
Unigene	335513
Immunogen	Recombinant fragment of human Factor XIIIa protein (aa46-181) (exact sequence is proprietary)
Antibody target cellular localization	Cytoplasmic, Secreted (extracellular)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, ELISA: 2-4 ug/mL for coating order Ab without BSA, Immunofluorescence: 0.5-1 ug/mL, Immunohistology (formalin) 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 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Western blotting 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user
Positive control	293T Cells. Placenta.
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
Cell/tissue expression	Platelets, Uterus, Megakaryocytes, Monocytes/macrophages, Placenta, Prostate
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.

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 25, 2024 at 07:28:07 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mouse-monoclonal-anti-factor-xiii-a-f13a11448/>