

FLI1 Monoclonal Mouse Antibody (FLI1/1312)

Product Description

This antibody recognizes a protein of 51 kDa, which is identified as FLI1. This protein, a member of the ETS family of DNA binding transcription factors, is involved in cellular proliferation and tumorigenesis. Ets-1 is the prototype member of a family of genes identified on the basis of homology to the v-Ets oncogene isolated from the E26 erythroblastosis virus. Members of the Ets gene family share a highly conserved carboxy-terminal domain containing a sequence related to the SV40 large T antigen nuclear localization signal sequence. Approximately 90% of Ewing s Sarcoma (EWS)/Primitive Neuroectodermal Tumors (PNET) have a specific translocation, t(11;22)(q24;q12), which results in fusion of EWS to Fli-1, and production of an EWS-Fli-1 fusion protein. Among normal tissues only endothelial cells and small lymphocytes express Fli-1. This protein is expressed in majority of vascular tumors including angiosarcomas, hemangioendotheliomas, hemangiomas, and Kaposi's Sarcomas. High sensitivity and specificity of Fli-1 equals to or exceeds that of the established vascular markers like CD31, CD34, and Factor VIII.

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 1312, Anti-FLI1 (FLI1/1312)

Product attributes

Antibody number	#1312
Antibody reactivity (target)	FLI1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	FLI1/1312
Isotype	IgG
Molecular weight	51 kDa
Synonyms	ERGB transcription factor; Ewing Sarcoma breakpoint region 2 (EWSR2); FLI1; FLI1 EWS fusion gene; Friend leukemia integration 1 (FLI1) transcription factor; Friend leukemia virus integration 1; Proto-oncogene Fli-1; SIC1; Transcription factor ERGB; Viral integration region FLI1
Human gene symbol	FLI1
Entrez gene ID	2313
SwissProt	Q01543
Unigene	504281
Immunogen	Recombinant fragment (aa 107-261) of human FLI1 protein (exact sequence is proprietary)
Verified antibody applications	WB (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Antibody application notes	Western blotting 1-2 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Immunofluorescence 1-2 ug/mL, Optimal dilution for a specific application should be determined by user
Positive control	THP1 and Raji cells. Tonsil, Ewing s sarcoma (EWS) and Angiosarcoma
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
Antibody research areas	Cancer, Immunology, Transcription factors
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.
Cell/tissue expression	Endothelial cells
Tumor expression	Sarcoma

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 July 14, 2025 at 01:03:35 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mouse-monoclonal-anti-fli1-fli11312/>