

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
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
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, LI-COR Bioscience.