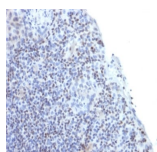


ETS1 Monoclonal Mouse Antibody (ETS1/1282)



Product Description

ETS1 proto-oncogene is an important transcription factor that plays a role in cell proliferation and differentiation. ETS1 is related to the growth of carcinoma cells by its regulation of the transcription of matrix metalloproteinases and urokinase-type plasminogen activator. The processes of tumor invasion and metastasis depend on the increased proteolytic activity of the invading tumor cells that may involve matrix metalloproteinases, cathepsins B and D and plasminogen activator in the metastatic cascade. ETS1 is preferentially expressed in lymphoid cells, where it is essential for the maintenance of the normal pool of resting T and B cells. ETS1 expression level and distribution are differentially controlled in resting, activated and apoptotic lymphocytes.

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 1282, Anti-ETS1- (ETS1/1282)

Product attributes

Antibody number	#1282
Antibody reactivity (target)	ETS1-
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ETS1/1282
Isotype	IgG
Molecular weight of antigen	54 kDa
Synonyms	Avian erythroblastosis virus E26 (v ets) oncogene homolog 1; ETS1; EWSR2; p54; v-ets erythroblastosis virus E26 oncogene homolog 1
Human gene symbol	ETS1
Entrez gene ID	2113
SwissProt	P14921
Unigene	369438
Immunogen	Recombinant human ETS1 fragment (aa137-230) (exact sequence is proprietary)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	HepG2 and HeLa cells. Tonsil, Skin, Pancreas, Kidney and Lymphoma.
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, 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.

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 June 13, 2026 at 06:19:32 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/mouse-monoclonal-anti-ets1-ets11282/>