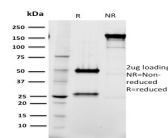


# Tal1 Monoclonal Mouse Antibody (BTL73)



## Product Description

Activation of TAL1 characterizes up to 60% of cases of human T cell acute lymphoblastic leukemia, making it the most frequent gain-of-function mutation observed in this disorder. TAL1 (also designated SCL) is a serine phosphoprotein and basic helix-loop-helix transcription factor known to regulate embryonic hematopoiesis. This transcription factor binds as a heterodimer with E2A and HEB/HTF4 to a nucleotide sequence motif termed the E-box. In addition, leukemogenesis is accelerated dramatically by transgenic co-expression of TAL1 and the catalytic subunit of casein kinase IIHLH transcription factors.

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 2852, Anti-Tal1 (BTL73)**

## Product attributes

<b>Antibody number</b>	#2852
<b>Antibody reactivity (target)</b>	Tal1
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	BTL73
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight</b>	42 kDa
<b>Synonyms</b>	bHLHa17; Class A basic helix-loop-helix protein 17; OTTHUMP0000009563; OTTHUMP0000009564; SCL; STEM CELL LEUKEMIA HEMATOPOIETIC TRANSCRIPTION FACTOR; Stem cell protein; T cell acute lymphocytic leukemia 1; T cell acute lymphocytic leukemia 1 protein; T cell acute lymphocytic leukemia 1 protein; T cell leukemia/lymphoma 5 protein; T-cell acute lymphocytic leukemia protein 1; T-cell leukemia/lymphoma protein 5; Tal 1 product; TAL 1 protein; TAL bHLH transcription factor 1 erythroid differentiation factor; TAL-1
<b>Human gene symbol</b>	TAL1
<b>Entrez gene ID</b>	6886
<b>SwissProt</b>	P17542
<b>Unigene</b>	705618
<b>Immunogen</b>	Recombinant human full-length TAL1 protein
<b>Antibody target cellular localization</b>	Nucleus
<b>Species reactivity</b>	Human
<b>Expected antibody applications</b>	IHC (published for clone), ChIP (published for clone), EMSA (published for clone), Flow (intracellular) (published for clone), IHC (FFPE) (published for clone)
<b>Antibody application notes</b>	ELISA: For coating, order antibody without BSA; Western blot 0.5-1.0 ug/mL; 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
<b>Positive control</b>	A549 xenograft.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	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	<a href="#">CF@405S Features</a>
BNC88	CF@488A	490/515	488	GFP, FITC	<a href="#">CF@488A Features</a>
BNC68	CF@568	562/583	532, 561	RFP, TRITC	<a href="#">CF@568 Features</a>
BNC94	CF@594	593/614	561	Texas Red®	<a href="#">CF@594 Features</a>
BNC40	CF@640R	642/662	633-640	Cy@5	<a href="#">CF@640R Features</a>
BNC47	CF@647	650/665	633-640	Cy@5	<a href="#">CF@647 Features</a>
BNC74	CF@740	742/767	633-685	775/50	<a href="#">CF@740 Features</a>
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. EMBO J (1997) 16:3145-3157. (WB)
2. Br J Haematol (1998) 102: 449-457. (Flow; IHC)
3. Blood (2002) 100(2): 491-500. (WB; IHC, FFPE; EMSA)
4. Oncogene (2011) 30: 1252-1260. (ChIP)
- 5.

This datasheet was generated on April 24, 2025 at 12:02:51 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/tal1-monoclonal-mouse-antibody-btl73/>