Spermidine or Spermine N1-Acetyltransferase 1 Monoclonal Mouse Antibody (CPTC-SAT1-3)



Product Description

Spermidine/spermine N1-acetyltransferase 1 (SAT1 or SSAT1) is the key regulatory enzyme in the catabolism of polyamines, catalyzing acetylation of spermidine or spermine to generate N1-acetyl spermidine or N1-acetyl spermine, and N1, N12-diacetylspermine. The cellular level of SAT1 is normally extremely low, but it is induced rapidly by a variety of stimuli, including polyamines, polyamine analogs, toxic chemicals, certain drugs, and growth factors. Downregulation of SAT1 has been reported in Epstein-Barr virus positive Burkitt's lymphoma cells.

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 2273, Anti-Spermidine or Spermine N1-Acetyltransferase 1 (CPTC-SAT1-3)

Product attributes

Product attributes				
Antibody number	1 2273			
Antibody reactivity (target)	imary			
Antibody type	Primary			
Host species	permidine or Spermine N1-Acetyltransferase 1 firmary pouse procolonal pTC-SAT1-3 G2a, kappa D kDa C21; Diamine acetyltransferase 1; Diamine N acetyltransferase EC 2.3.1.57; KDSD; KSDX; Polyamine N acetyltransferase 1; utrescine acetyltransferase; SAT1; spermidine/spermine N1 petyltransferase; SAT1; spermidine/spermine N1 petyltransferase alpha; SSAT; SSAT1 AT1 AT1 AT1 BAT1 BAT1 BAT1 BAT1 BAT1			
Clonality	Monoclonal			
Clone	CPTC-SAT1-3			
Isotype	IgG2a, kappa			
Molecular weight	20 kDa			
Synonyms	DC21; Diamine acetyltransferase 1; Diamine N acetyltransferase 1; EC 2.3.1.57; KDSD; KSDX; Polyamine N acetyltransferase 1; Putrescine acetyltransferase; SAT1; spermidine/spermine N1 acetyltransferase alpha; SSAT; SSAT1			
Human gene symbol	SAT1			
Entrez gene ID	6303			
SwissProt	P21673			
Unigene	28491			
Immunogen	Recombinant full-length human SAT1 protein			
Antibody target cellular localization	Cytoplasmic			
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	Human liver tissue lysate; human tonsil or kidney.			
Shipping condition	Room temperature			
Storage Conditions	Store at 2 to 8 $^{\circ}$ C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ C			
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	PBS without azide Monospecific			
Validated in protein array Shelf life				
, ,	Monospecific Guaranteed for at least 24 months from date of receipt when			

Call us: 800-304-5357 Email: btinfo@biotium.com

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 4, 2025 at 06:41:45 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/spermidine-or-spermine-n1-acetyltransferase-1-monoclonal-mouse-antibody-cptc-sat1-3/