

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

| kDa | | NF | | R |
|------|---|----|---|--------------------|
| 250 | | | | |
| 150 | | - | | |
| 100 | | | | |
| 75 — | - | | | 2ug loadir |
| 50 | - | | - | NR=Non- reduced |
| 37 — | - | | | R=reduced |
| 25 | - | | _ | |
| 20 | | | | |
| 15 | - | | | |

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)

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

| Product attributes | | | |
|--|---|--|--|
| Antibody number | #2273 | | |
| Antibody reactivity (target) | Spermidine or Spermine N1-Acetyltransferase 1 | | |
| Antibody type | Primary | | |
| Host species | Mouse | | |
| Clonality | Monoclonal | | |
| Clone | CPTC-SAT1-3 | | |
| Isotype | lgG2a, kappa | | |
| Molecular weight | 20 kDa | | |
| Synonyms | DC21; Diamine acetyltransferase 1; Diamine N acetyltransferase 1; EC 2.3.1.57; KOSD; 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 | | |
| Species reactivity Antibody application notes | Human For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody, oonjugates 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 | | |
| | 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 | | |
| Antibody application notes | For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody, opiupaties 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 | | |
| Antibody application notes Positive control | For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody, oonjugates 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 Human liver tissue lysate; human tonsil or kidney. | | |
| Antibody application notes Positive control Shipping condition | For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary 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 Human liver tissue lysate; human tonsil or kidney. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, | | |
| Antibody application notes Positive control Shipping condition Storage Conditions | For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody, Ooptimal 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 Human liver tissue lysate; human tonsil or kidney. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C | | |
| Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate | For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary 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 Human liver tissue lysate; human tonsil or kidney. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA/0.05% azide, HRP | | |
| Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation | For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary 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 Human liver tissue lysate; human tonsil or kidney. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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 application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Validated in protein array | For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary 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 Human liver tissue lysate; human tonsil or kidney. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP 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 Monospecific Guaranteed for at least 24 months from date of receipt when | | |

| 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 12, 2025 at 02:36:35 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/