

PRODUCT INFORMATION



8-Bu(3-yne)T-NAD⁺ (sodium salt)

Item No. 38477

CAS Registry No.: 2022926-15-2

Formal Name: 8-(3-butyn-1-ylthio)-adenosine 5'-(trihydrogen diphosphate), P'→5'-ester with 3-(aminocarbonyl)-1-β-D-ribofuranosylpyridinium, inner salt, compd. with N,N-diethylethanamine, monosodium salt

Synonyms: Click Tag™ 8-Bu(3-yne)T-NAD, 8-Bu(3-yne)T-Nicotinamide adenine dinucleotide, β-Nicotinamide-8-(3-butynylthio)adenine dinucleotide

MF: C₂₅H₃₀N₇O₁₄P₂S • Na

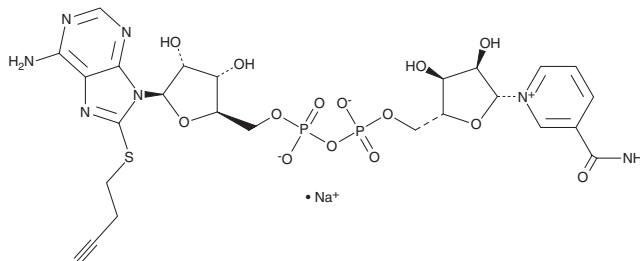
FW: 769.5

Purity: ≥95%

Supplied as: A solid

Storage: -80°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

8-Bu(3-yne)T-NAD⁺ (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 8-Bu(3-yne)T-NAD⁺ (sodium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

8-Bu(3-yne)T-NAD⁺ is a clickable form of the signaling molecule and enzyme cofactor NAD⁺ (Item No. 16077).^{1,2} It has been used in the determination of poly(ADP-ribose) polymerase 1- (PARP1-), PARP2-, and PARP3-dependent genomic ADP-ribosylation sites and PARP7 protein substrates using analog-sensitive gatekeeper mutants of these proteins.

References

1. Gibson, B.A., Zhang, Y., Jiang, H., *et al.* Chemical genetic discovery of PARP targets reveals a role for PARP-1 in transcription elongation. *Science* **353**(6294), 45-50 (2016).
2. Parsons, L.H.P., Challa, S., Gibson, B.A., *et al.* Identification of PARP-7 substrates reveals a role for MARYlation in microtubule control in ovarian cancer cells. *Elife* **10**, e60481 (2021).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM