

PRODUCT INFORMATION



8-bromo NAD⁺ (sodium salt)

Item No. 38475

CAS Registry No.: 2022926-16-3

Formal Name: 8-bromo-adenosine 5'-(trihydrogen diphosphate), P'→5'-ester with 3-(aminocarbonyl)-1-β-D-ribofuranosylpyridinium, inner salt, monosodium salt

Synonyms: N(8-bromo-A)D⁺, 8-bromo Nicotinamide adenine dinucleotide

MF: C₂₁H₂₅BrN₇O₁₄P₂ • Na

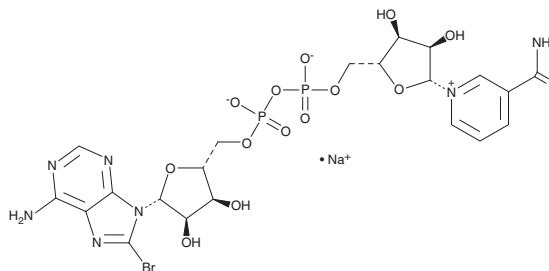
FW: 764.3

Purity: ≥98%

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

Description

8-bromo NAD⁺ is a prodrug form of the cyclic ADP-ribose (cADPR) inhibitor 8-bromo cADPR (Item No. 18442).¹ It is converted to 8-bromo-cADPR by CD38. 8-bromo NAD⁺ (1 mM) prevents increases in intracellular calcium levels and chemotaxis induced by N-formyl-Met-Leu-Phe (fMLP; Item No. 21495) in isolated mouse bone marrow-derived neutrophils. It also decreases LPS-induced nitrite production and TNF-α and IL-2 secretion in mouse primary microglial cells when used at a concentration of 100 μM.²

References

1. Partida-Sánchez, S., Cockayne, D.A., Monard, S., *et al.* Cyclic ADP-ribose production by CD38 regulates intracellular calcium release, extracellular calcium influx and chemotaxis in neutrophils and is required for bacterial clearance *in vivo*. *Nat. Med.* **7**(11), 1209-1216 (2001).
2. Mayo, L., Jacob-Hirsch, J., Amariglio, N., *et al.* Dual role of CD38 in microglial activation and activation-induced cell death. *J. Immunol.* **181**(1), 92-103 (2008).

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.

Copyright Cayman Chemical Company, 04/17/2026

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