

PRODUCT INFORMATION



7-deaza-8-bromo NAD⁺ (sodium salt)

Item No. 38479

Formal Name: 3-(aminocarbonyl)-1-[5-O-[hydroxy(phosphonooxy)phosphinyl]-β-D-ribofuranosyl]-pyridinium, P'→5'-ester with 6-bromo-7-β-D-ribofuranosyl-7H-pyrrolo[2,3-d]pyrimidin-4-amine, inner salt, monosodium salt

Synonym: 7-deaza-8-bromo Nicotinamide adenine dinucleotide

MF: C₂₂H₂₆BrN₆O₁₄P₂ • Na

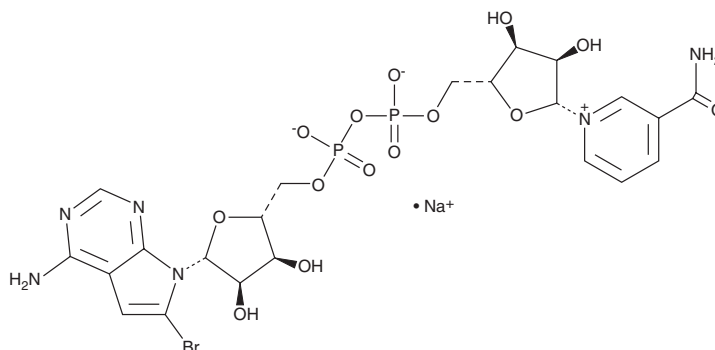
FW: 763.3

Purity: ≥95%

Supplied as: A solid

Storage: -80°C

Stability: ≥2 years



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

Description

7-deaza-8-bromo NAD⁺ is a brominated and deazaadenine derivative of the signaling molecule and cofactor NAD⁺ (Item No. 16077).¹ It has been used as a synthetic intermediate in the synthesis of cyclic and membrane-permeable diphosphate modulators of transient receptor potential melastatin 2 (TRPM2).

Reference

1. Moreau, C., Kirchberger, T., Swarbrick, J.M., *et al.* Structure-activity relationship of adenosine 5'-diphosphoribose at the transient receptor potential melastatin 2 (TRPM2) channel: Rational design of antagonists. *J. Med. Chem.* **56**(24), 10079-100102 (2013).

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, 05/24/2023

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