

# PRODUCT INFORMATION



## 8-cPeT NAD<sup>+</sup> (sodium salt)

Item No. 38480

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

**Synonym:** 8-cPeT Nicotinamide adenine dinucleotide

**MF:** C<sub>26</sub>H<sub>34</sub>N<sub>7</sub>O<sub>14</sub>P<sub>2</sub>S • Na

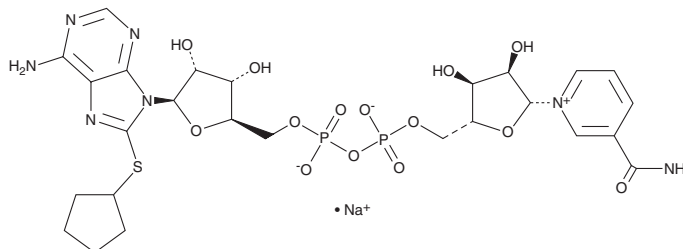
**FW:** 785.6

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

### Description

8-cPeT NAD<sup>+</sup> is a derivative of the signaling molecule and enzyme cofactor NAD<sup>+</sup> (Item No. 16077).<sup>1</sup> It has been used in the screening of analog-sensitive poly(ADP-ribose) polymerase 1 (PARP1) gatekeeper mutations.

### Reference

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).

#### 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, 04/20/2026

#### 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](http://WWW.CAYMANCHEM.COM)