

# PRODUCT INFORMATION



## 7-Deazaadenosine-5'-O-diphosphate (sodium salt)

Item No. 38451

**Formal Name:** ((2R,3S,4R,5R)-5-(4-amino-7H-pyrrolo[2,3-d]pyrimidin-7-yl)-3,4-dihydroxytetrahydrofuran-2-yl)methyl diphosphate, trisodium salt

**Synonyms:** 7-CH-ADP, Tubercidin 5'-diphosphate, TuDP

**MF:** C<sub>11</sub>H<sub>13</sub>N<sub>4</sub>O<sub>10</sub>P<sub>2</sub> • 3Na

**FW:** 492.2

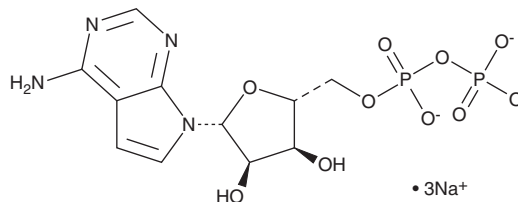
**Purity:** ≥95%

**Supplied as:** A solution in water

**Storage:** -80°C

**Stability:** ≥2 years

**Item Origin:** Synthetic



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

### Description

7-Deazaadenosine-5'-O-diphosphate is a modified nucleotide and metabolite of tubercidin (Item No. 35834).<sup>1</sup> It is formed from tubercidin *via* phosphorylation, and is further phosphorylated to the active metabolite 7-deazaadenosine-5'-O-triphosphate (Item No 38377).

### Reference

1. Ross, A.F. and Jaffe, J.J. Effects of tubercidin and its ribonucleotides on various metabolic pathways in *Schistosoma mansoni*. *Biochem. Pharmacol.* **21(22)**, 3059-3069 (1972).

#### 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, 05/05/2023

#### 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