

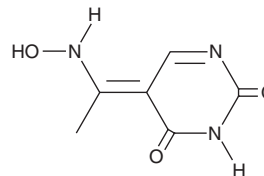
# PRODUCT INFORMATION



**NSC 232003**

Item No. 38523

**CAS Registry No.:** 1905453-18-0  
**Formal Name:** 5-[1-(hydroxyamino)ethylidene]-2,4(3H,5H)-pyrimidinedione  
**MF:** C<sub>6</sub>H<sub>7</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 169.1  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

## Laboratory Procedures

NSC 232003 is supplied as a solid. A stock solution may be made by dissolving the NSC 232003 in water. We do not recommend storing the aqueous solution for more than one day.

## Description

NSC 232003 is an inhibitor of the protein-protein interaction between E3 ubiquitin-protein ligase UHRF1 and DNA methyltransferase 1 (DNMT1; IC<sub>50</sub> = 15 μM in U251 glioma cells).<sup>1</sup> It decreases 5-methylcytosine (5-mC) levels in U251 cells (IC<sub>50</sub> = 15 μM).

## Reference

1. Myrianthopoulos, V., Cartron, P.F., Liutkevičiūtė, Z., *et al.* Tandem virtual screening targeting the SRA domain of UHRF1 identifies a novel chemical tool modulating DNA methylation. *Eur. J. Med. Chem.* **114**, 390-396 (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, 05/04/2023

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