

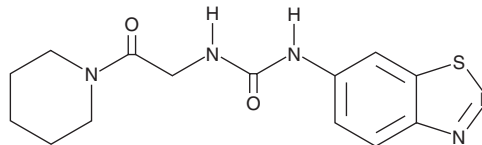
# PRODUCT INFORMATION



## UNC2327

Item No. 18122

**CAS Registry No.:** 1426152-53-5  
**Formal Name:** N-1,2,3-benzothiadiazol-6-yl-N'-[2-oxo-2-(1-piperidinyl)ethyl]-urea  
**MF:** C<sub>14</sub>H<sub>17</sub>N<sub>5</sub>O<sub>2</sub>S  
**FW:** 319.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 256, 311 nm  
**Supplied as:** A solid  
**Storage:** Room temperature  
**Stability:** ≥4 years



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

### Laboratory Procedures

UNC2327 is supplied as a solid. A stock solution may be made by dissolving the UNC2327 in the solvent of choice, which should be purged with an inert gas. UNC2327 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of UNC2327 in these solvents is approximately 2 and 10 mg/ml, respectively.

### Description

UNC2327 is an allosteric inhibitor of protein arginine methyltransferase 3 (PRMT3; IC<sub>50</sub> = 230 nM).<sup>1</sup>

### Reference

1. Liu, F., Li, F., Ma, A., *et al.* Exploiting an allosteric binding site of PRMT3 yields potent and selective inhibitors. *J. Med. Chem.* **56**(5), 2110-2124 (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, 10/05/2022

#### 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