

# PRODUCT INFORMATION



## Cdk9 Inhibitor

Item No. 42512

**CAS Registry No.:** 1415559-43-1  
**Formal Name:** N-[5-[[6-[3-(1,3-dihydro-1,3-dioxo-2H-isoindol-2-yl)phenyl]-4-pyrimidinyl]amino]-2-methylphenyl]-methanesulfonamide

**Synonyms:** Cdk9-IN-1, Cyclin-dependent Kinase 9 Inhibitor, Cyclin-dependent Kinase 9 Inhibitor-1, Cyclin-dependent Kinase 9-IN-1, Cdk9 Inhibitor 1, KIN001-270

**MF:** C<sub>26</sub>H<sub>21</sub>N<sub>5</sub>O<sub>4</sub>S

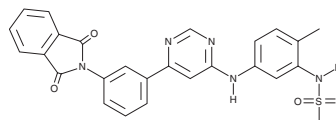
**FW:** 499.5

**Purity:** ≥95%

**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

Cdk9 inhibitor is supplied as a solid. A stock solution may be made by dissolving the Cdk9 inhibitor in the solvent of choice, which should be purged with an inert gas. Cdk9 inhibitor is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

Cdk9 inhibitor is an inhibitor of cyclin-dependent kinase 9 (Cdk9; IC<sub>50</sub> = 39 nM).<sup>1</sup> It is selective for Cdk9 over Cdk1-7 (IC<sub>50</sub>s = >1,000 nM for all). Cdk9 inhibitor (6-200 nM) inhibits porcine reproductive and respiratory syndrome virus (PRRSV) replication in PRRSV-infected primary porcine alveolar macrophages.<sup>2</sup>

### References

1. Németh, G., Varga, Z., Greff, Z., *et al.* Novel, selective CDK9 inhibitors for the treatment of HIV infection. *Curr. Med. Chem.* **18(3)**, 342-358 (2011).
2. Wang, M.D., Yang, L., Meng, J.J., *et al.* Functionally active cyclin-dependent kinase 9 is essential for porcine reproductive and respiratory syndrome virus subgenomic RNA synthesis. *Mol. Immunol.* **135**, 351-364 (2021).

#### 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, 12/19/2024

#### 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