

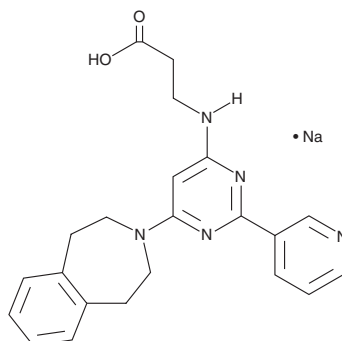
# PRODUCT INFORMATION



## GSK-J2 (sodium salt)

Item No. 12056

CAS Registry No.: 2108665-15-0  
Formal Name: 3-((6-(4,5-dihydro-1H-benzo[d]azepin-3(2H)-yl)-2-(pyridin-3-yl)pyrimidin-4-yl)amino)propanoic acid, monosodium salt  
MF:  $C_{22}H_{23}N_5O_2 \cdot Na$   
FW: 412.4  
Purity:  $\geq 90\%$   
UV/Vis.:  $\lambda_{max}$ : 242 nm  
Supplied as: A crystalline solid  
Storage:  $-20^\circ C$   
Stability:  $\geq 4$  years



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

### Description

The histone H3 lysine 27 (H3K27) demethylase JMJD3 plays important roles in the transcriptional regulation of cell differentiation, development, the inflammatory response, and cancer.<sup>1,2</sup> GSK-J2 is a pyridine regio-isomer of GSK-J1 which poorly inhibits JMJD3 ( $IC_{50} > 100 \mu M$ ), making it an appropriate negative control for *in vitro* studies involving GSK-J1.<sup>3</sup> For *in vivo* research, cell-permeable prodrug forms of GSK-J1 and GSK-J2 are available as GSK-J4 (Item No. 12073) and GSK-J5 (Item No. 12074), respectively. See the Structural Genomics Consortium (SGC) website for more information.

### References

1. Agger, K., Cloos, P.A.C., Christensen, J., *et al.* UTX and JMJD3 are histone H3K27 demethylases involved in HOX gene regulation and development. *Nature* **449**(7163), 731-4 (2011).
2. Hübner, M.R. and Spector, D.L. Role of H3K27 demethylases Jmjd3 and UTX in transcriptional regulation. Cold Spring Harb. Symp. Quant. Biol. **75**, 43-49 (2010).
3. Kruidenier, L., Chung, C.W., Cheng, Z., *et al.* A selective jumonji H3K27 demethylase inhibitor modulates the proinflammatory macrophage response. *Nature* **488**(7411), 404-408 (2012).

#### 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, 02/27/2024

#### 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