

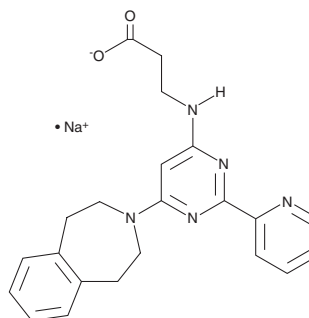
PRODUCT INFORMATION



GSK-J1 (sodium salt)

Item No. 12054

CAS Registry No.: 1797832-71-3
Formal Name: N-[2-(2-pyridinyl)-6-(1,2,4,5-tetrahydro-3H-3-benzazepin-3-yl)-4-pyrimidinyl]-β-alanine, monosodium salt
MF: C₂₂H₂₂N₅O₂ • Na
FW: 411.4
Purity: ≥95%
UV/Vis.: λ_{max}: 253 nm
Supplied as: A crystalline 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

GSK-J1 (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK-J1 (sodium salt) in the solvent of choice, which should be purged with an inert gas. GSK-J1 (sodium salt) is slightly soluble in DMSO.

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.^{1,2} GSK-J1 is a potent inhibitor of the H3K27 histone demethylases JMJD3 and UTX (IC₅₀s = 18 and 56 μM, respectively as measured by mass spectrometry; IC₅₀ = 60 nM in JMJD3 antibody-based assays).³ It is inactive against a panel of additional JMJD family demethylases, including several variants of JMJD2 and JMJD1 and, at concentrations up to 30 μM, has no effect on more than 100 different kinases or other unrelated proteins, including other chromatin-modifying enzymes such as histone deacetylases.³ Since the highly polar carboxylate group of GSK-J1 restricts its cellular permeability, a prodrug ethyl ester, GSK-J4 (Item No. 12073) has also been developed. 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. *Nat. Lett.* **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**, (2011).
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