

PRODUCT INFORMATION



GSK591

Item No. 18354

CAS Registry No.: 1616391-87-7

Formal Name: 2-(cyclobutylamino)-N-[(2S)-3-(3,4-dihydro-2(1H)-isoquinoliny)-2-hydroxypropyl]-4-pyridinecarboxamide

Synonyms: EPZ015866, GSK3203591

MF: $C_{22}H_{28}N_4O_2$

FW: 380.5

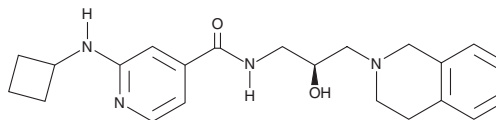
Purity: $\geq 95\%$

UV/Vis.: λ_{max} : 251, 340 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

GSK591 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK591 in the solvent of choice, which should be purged with an inert gas. GSK591 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GSK591 in ethanol is approximately 12 mg/ml and approximately 16 mg/ml in DMSO and DMF.

GSK591 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK591 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK591 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GSK591 is a chemical probe for protein arginine methyltransferase 5 (PRMT5) that potently inhibits the PRMT5/MEP50 complex from methylating histone H4 ($IC_{50} = 11$ nM) *in vitro*. GSK591 is selective for PRMT5 (up to 50 μM) relative to a panel of other methyltransferases. In Z-138 cells, GSK591 inhibits the symmetric arginine methylation of SmD3 with an EC_{50} value of 56 nM. For more information on GSK591 please visit the Structural Genomics Consortium (SGC). The negative control, SGC2096, for GSK591 is also available exclusively through the SGC.

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, 12/06/2022

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