

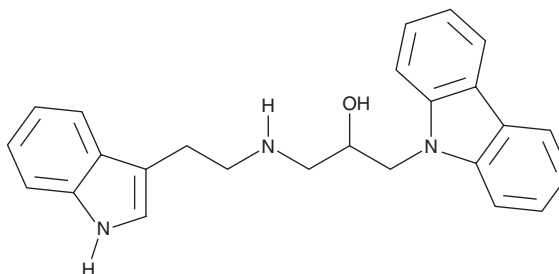
PRODUCT INFORMATION



DC-05

Item No. 36584

CAS Registry No.: 890643-16-0
Formal Name: α -[[[2-(1H-indol-3-yl)ethyl]amino]methyl]-9H-carbazole-9-ethanol
MF: $C_{25}H_{25}N_3O$
FW: 383.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 223, 235, 261, 292 nm
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

DC-05 is supplied as a solid. A stock solution may be made by dissolving the DC-05 in the solvent of choice, which should be purged with an inert gas. DC-05 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of DC-05 in these solvents is approximately 11 and 12 mg/ml, respectively. DC-05 is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of DC-05 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of DC-05 in PBS (pH 7.2) is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

DC-05 is an inhibitor of DNA methyltransferase 1 (DNMT1; $IC_{50} = 10.3 \mu\text{M}$).¹ It is selective for DNMT1 over DNMT3A and DNMT3B (IC_{50} s = $>200 \mu\text{M}$), as well as the additional methyltransferases G9a, SUV39H1, MLL1, SET7/9, and PRMT1 (IC_{50} s = >150 , >150 , >150 , >150 , and $37.1 \mu\text{M}$, respectively). DC-05 (2.5 - $10 \mu\text{M}$) reduces proliferation of HCT116 colon cancer cells.

Reference

1. Chen, S., Wang, Y., Zhou, W., *et al.* Identifying novel selective non-nucleoside DNA methyltransferase 1 inhibitors through docking-based virtual screening. *J. Med. Chem.* **57**(21), 9028-9041 (2014).

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, 12/20/2022

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