

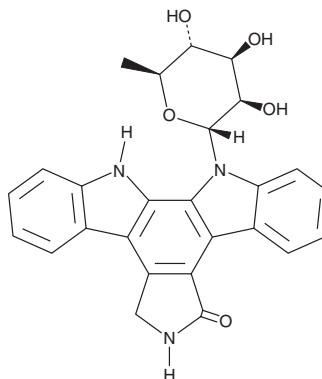
PRODUCT INFORMATION



K252d

Item No. 28469

CAS Registry No.: 105114-22-5
Formal Name: 13-(6-deoxy- α -L-mannopyranosyl)-
6,7,12,13-tetrahydro-5H-indolo[2,3-a]
pyrrolo[3,4-c]carbazol-5-one
MF: $C_{26}H_{23}N_3O_5$
FW: 457.5
Purity: $\geq 70\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Bacterium/*Nonomuraea longicatena*



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

Laboratory Procedures

K252d is supplied as a solid. A stock solution may be made by dissolving the K252d in the solvent of choice, which should be purged with an inert gas. K252d is soluble in organic solvents such as ethanol and DMSO.

Description

K252d is an indolocarbazole alkaloid found in *Nocardiosis*.¹ It is a PKC inhibitor that inhibits PKC isolated from rat brain ($IC_{50} = 350$ nM). It also inhibits calcium- and calmodulin-dependent phosphodiesterase isolated from bovine heart ($IC_{50} = 46.2$ μM).

Reference

1. Nakanishi, S., Matsuda, Y., Iwahashi, K., et al. K-252b, c and d, potent inhibitors of protein kinase C from microbial origin. *J. Antibiot. (Tokyo)* **39**(8), 1066-1071 (1986).

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, 08/16/2023

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