

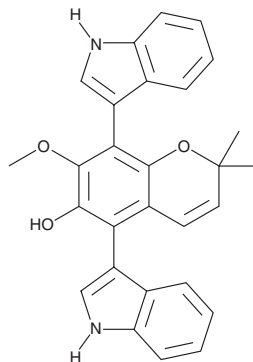
PRODUCT INFORMATION



Kumbicin C

Item No. 25015

CAS Registry No.: 1878151-58-6
Formal Name: 5,8-di-1H-indol-3-yl-7-methoxy-2,2-dimethyl-2H-1-benzopyran-6-ol
MF: $C_{28}H_{24}N_2O_3$
FW: 436.5
Purity: $\geq 95\%$
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

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

Description

Kumbicin C is a *bis*-indolyl benzenoid fungal metabolite produced by *A. kumbius* FRR6049.¹ It inhibits the growth of NS-1 mouse myeloma cells ($IC_{50} = 0.74 \mu\text{g/ml}$). Kumbicin C also inhibits the growth of the Gram-positive bacteria *B. subtilis* (MIC = $1.6 \mu\text{g/ml}$).

Reference

1. Lacey, H.J., Vuong, D., Pitt, J.I., *et al.* Kumbicins A–D: Bis-Indolyl benzenoids and benzoquinones from an Australian soil fungus, *Aspergillus kumbius*. *Aust. J. Chem.* **69**(2), 152-160 (2015).

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, 12/19/2022

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