

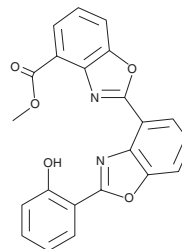
PRODUCT INFORMATION



UK 1

Item No. 20139

CAS Registry No.: 151271-53-3
Formal Name: 2'-(2-hydroxyphenyl)-[2,4'-bibenzoxazole]-4-carboxylic acid, methyl ester
MF: C₂₂H₁₄N₂O₅
FW: 386.4
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

UK 1 is supplied as a crystalline solid. A stock solution may be made by dissolving the UK 1 in the solvent of choice, which should be purged with an inert gas. UK 1 is soluble in organic solvents such as DMSO and dimethyl formamide.

Description

UK 1 is a metabolite from *Streptomyces* sp. 517-02 that demonstrates cytotoxic activity against various cancer cell lines, though it does not have antibacterial effects.¹ It has been shown to inhibit topoisomerase II and also to inhibit hepatitis C viral replication.²⁻³

References

1. Ueki, M., Ueno, K., Miyadoh, S., *et al.* UK-1, a novel cytotoxic metabolite from *Streptomyces* sp. 517-02. I. Taxonomy, fermentation, isolation, physico-chemical and biological properties. *J. Antibiot (Tokyo)*. **46(7)**, 1089-1094 (1993).
2. Reynolds, M. B., DeLuca, M. R., and Kerwin, S. M. The novel Bis(benzoxazole) Cytotoxic natural product UK-1 is a magnesium ion-dependent DNA binding agent and inhibitor of human topoisomerase II. *Bioorg. Chem.* **27(4)**, 326-337 (1999).
3. Ward, D. N., Talley, D. C., Tavag, M., *et al.* UK-1 and structural analogs are potent inhibitors of hepatitis C virus replication. *Bioorg. Med. Chem. Lett.* **24(2)**, 609-612 (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, 11/11/2022

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