

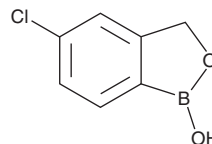
PRODUCT INFORMATION



AN2718

Item No. 20708

CAS Registry No.: 174672-06-1
Formal Name: 5-chloro-1,3-dihydro-1-hydroxy-2,1-benzoxaborole
MF: C₇H₆BClO₂
FW: 168.4
Purity: ≥98%
UV/Vis.: λ_{max}: 228, 271, 279 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

AN2718 is supplied as a crystalline solid. A stock solution may be made by dissolving the AN2718 in the solvent of choice, which should be purged with an inert gas. AN2718 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AN2718 in these solvents is approximately 30 mg/ml.

AN2718 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AN2718 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. AN2718 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AN2718 is a benzoxaborole compound that displays broad antifungal effectiveness *in vitro*.¹ It inhibits leucyl-tRNA synthetase (LeuRS) from molds and yeasts (IC₅₀s = 2 and 4.2 μM for *A. fumigatus* and *C. albicans* LeuRS, respectively), blocking protein synthesis.¹

Reference

1. Seiradake, E., Ip, E., Rock, F., *et al.* Antifungal activity and mechanism of action of a benzoxaborole, AN2718, which is in development for the treatment of tinea pedis. *48th Annual ICAAC/46th Annual IDSA* (2008).

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, 12/16/2022

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