

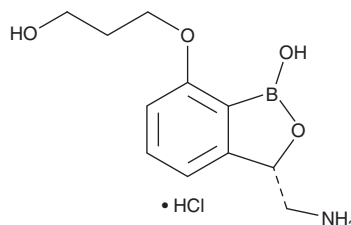
# PRODUCT INFORMATION



## AN3365 (hydrochloride)

Item No. 20572

**CAS Registry No.:** 1234563-16-6  
**Formal Name:** 3-[[[(3S)-3-(aminomethyl)-1,3-dihydro-1-hydroxy-2,1-benzoxaborol-7-yl]oxy]-1-propanol, monohydrochloride  
**Synonyms:** Epetraborole, GSK2251052  
**MF:** C<sub>11</sub>H<sub>16</sub>BNO<sub>4</sub> • HCl  
**FW:** 273.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 230, 284 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

AN3365 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the AN3365 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. AN3365 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

### Description

AN3365 is an inhibitor of leucyl-tRNA synthetase (LeuRS; IC<sub>50</sub> = 0.31 μM for the *E. coli* enzyme).<sup>1</sup> It is active against a variety of Gram-negative and Gram-positive bacteria, including *E. coli*, *E. faecalis*, *S. epidermidis*, *S. pneumoniae*, and *S. pyogenes* (MICs = 4, 8, 2, 2, and 2 μg/ml, respectively), as well as a panel of multidrug-resistant (MDR) clinical isolates of the Gram-negative bacteria *A. baumannii*, *Enterobacteriaceae*, and *P. aeruginosa* (MICs = 0.5-4, 0.5-2, and 1-4 μg/ml, respectively). AN3365 inhibits protein, RNA, DNA, and fatty acid synthesis in *E. coli* (IC<sub>50</sub>s = 8.3, 105, 366, and 396 μg/ml, respectively). It reduces the number of colony forming units (CFUs) in neutropenic mouse thigh models of *E. coli* or *P. aeruginosa* infection when administered at a dose of 30 mg/kg.

### Reference

1. Hernandez, V., Crépin, T., Palencia, A., *et al.* Discovery of a novel class of boron-based antibacterials with activity against gram-negative bacteria. *Antimicrob. Agents and Chemother.* **57**(3), 1394-1403 (2013).

#### 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, 10/18/2021

#### 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