

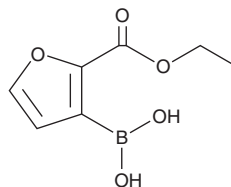
PRODUCT INFORMATION



TP-4748

Item No. 31459

CAS Registry No.: 1150114-62-7
Formal Name: 3-borono-2-furancarboxylic acid, 2-ethyl ester
MF: C₇H₉BO₅
FW: 184.0
Purity: ≥98%
UV/Vis.: λ_{max}: 258 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

TP-4748 is supplied as a crystalline solid. A stock solution may be made by dissolving the TP-4748 in the solvent of choice, which should be purged with an inert gas. TP-4748 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TP-4748 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

TP-4748 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TP-4748 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. TP-4748 has a solubility of approximately 0.50 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TP-4748 is a heterocyclic building block.¹ It has been used in the synthesis of prolyl-tRNA synthase inhibitors.

Reference

1. Lee, B.-Y., Cho, M.-J., Lee, H.-G., *et al.* Novel heterocyclic compound, its preparation method, and pharmaceutical composition comprising the same. *Deawoong Pharmaceutical Co., LTD.* **WO 2018/147626 A1** (2018).

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, 11/10/2022

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