

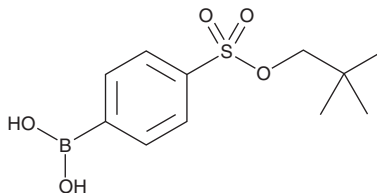
PRODUCT INFORMATION



4-(Neopentyloxysulfonyl)phenylboronic Acid

Item No. 21493

CAS Registry No.: 957060-74-1
Formal Name: 4-borono-benzenesulfonic acid,
1-(2,2-dimethylpropyl) ester
MF: C₁₁H₁₇BO₅S
FW: 272.1
Purity: ≥95%
UV/Vis.: λ_{max}: 229, 272 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

4-(Neopentyloxysulfonyl)phenylboronic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-(neopentyloxysulfonyl)phenylboronic acid in the solvent of choice, which should be purged with an inert gas. 4-(Neopentyloxysulfonyl)phenylboronic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 4-(neopentyloxysulfonyl)phenylboronic acid in these solvents is approximately 30 mg/ml.

4-(Neopentyloxysulfonyl)phenylboronic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-(neopentyloxysulfonyl)phenylboronic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 4-(Neopentyloxysulfonyl)phenylboronic acid has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

4-(Neopentyloxysulfonyl)phenylboronic acid is a monosulfonic acid that has been used as a building block in the synthesis of water-soluble dibenzothiophene (DBT) and dibenzothiophene S-oxide (DBTO) derivatives.¹

Reference

1. Omlid, S.M., Isor, A., Sulkowski, K.L., *et al.* Synthesis of aromatic disulfonic acids for water-soluble dibenzothiophene derivatives. *Synthesis* **50**(12), 2359-2366 (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, 11/11/2022

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