

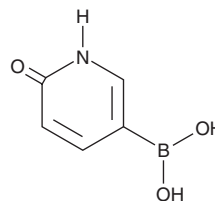
# PRODUCT INFORMATION



## 6-Hydroxypyridin-3-ylboronic Acid

Item No. 30492

**CAS Registry No.:** 903899-13-8  
**Formal Name:** B-(1,6-dihydro-6-oxo-3-pyridinyl)-boronic acid  
**MF:** C<sub>5</sub>H<sub>6</sub>BNO<sub>3</sub>  
**FW:** 138.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 243, 300 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

6-Hydroxypyridin-3-ylboronic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 6-hydroxypyridin-3-ylboronic acid in the solvent of choice, which should be purged with an inert gas. 6-Hydroxypyridin-3-ylboronic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 6-hydroxypyridin-3-ylboronic acid in these solvents is approximately 1 mg/ml in ethanol and 30 mg/ml in DMSO and DMF.

6-Hydroxypyridin-3-ylboronic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 6-hydroxypyridin-3-ylboronic acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 6-Hydroxypyridin-3-ylboronic acid has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

6-Hydroxypyridin-3-ylboronic acid is a heterocyclic building block.<sup>1,2</sup> It has been used in the synthesis of non-nucleoside inhibitors of hepatitis C virus (HCV) RNA-dependent RNA polymerase nonstructural protein 5B (NS5B).<sup>1</sup> 6-Hydroxypyridin-3-ylboronic acid has also been used in the synthesis of mammalian target of rapamycin (mTOR) inhibitors.<sup>2</sup>

### Reference

- Hendricks, R.T., Spencer, S.R., Blake, J.F., *et al.* 3-Hydroxyisoquinolines as inhibitors of HCV NS5B RNA-dependent RNA polymerase. *Bioorg. Med. Chem. Lett.* **19(2)**, 410-414 (2009).
- Verheijen, J.C., Richard, D.J., Curran, K., *et al.* Discovery of 4-morpholino-6-aryl-1H-pyrazolo[3,4-d]pyrimidines as highly potent and selective ATP-competitive inhibitors of the mammalian target of rapamycin (mTOR): Optimization of the 6-aryl substituent. *J. Med. Chem.* **52(24)**, 8010-8024 (2009).

#### 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, 05/05/2020

#### 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