

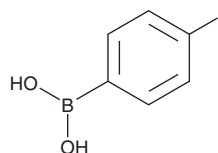
PRODUCT INFORMATION



(4-Iodophenyl)boronic Acid

Item No. 45679

CAS Registry No.: 5122-99-6
Formal Name: B-(4-iodophenyl)-boronic acid
Synonym: 4-IPBA
MF: $C_6H_6BIO_2$
FW: 247.8
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

(4-Iodophenyl)boronic acid is supplied as a solid. A stock solution may be made by dissolving the (4-iodophenyl)boronic acid in the solvent of choice, which should be purged with an inert gas. (4-Iodophenyl)boronic acid is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

(4-Iodophenyl)boronic acid is an enhancer of luminol (Item No. 16803) chemiluminescent peroxidation catalyzed by horseradish peroxidase (HRP).¹ HRP-induced chemiluminescent light emission increases by >100 -fold in the presence of (4-iodophenyl)boronic acid. (4-Iodophenyl)boronic acid has also been used as a fluorescent chemosensor in gold clusters for the detection of saccharides.²

References

1. Kricka, L.J., Cooper, M., and Ji, X. Synthesis and characterization of 4-iodophenylboronic acid: A new enhancer for the horseradish peroxidase-catalyzed chemiluminescent oxidation of luminol. *Anal. Biochem.* **240**, 119-125 (1996).
2. Thakarda, J., Dave, P., Bhowmik, S., et al. 4-Iodophenylboronic acid stabilized gold cluster as a new fluorescent chemosensor for saccharides based on excimer emission quenching. *J. Fluoresc.* **31(2)**, 441-454 (2021).

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, 06/01/2026

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