

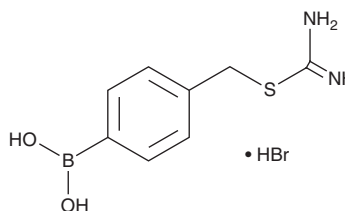
PRODUCT INFORMATION



BC-11 (hydrobromide)

Item No. 39522

CAS Registry No.: 443776-49-6
Formal Name: carbamimidothioic acid, (4-boronophenyl) methyl ester, monohydrobromide
Synonyms: uPA Inhibitor BC-11, Urokinase-type Plasminogen Activator Inhibitor BC-11
MF: $C_8H_{11}BN_2O_2S \cdot HBr$
FW: 291.0
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

BC-11 (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the BC-11 (hydrobromide) in the solvent of choice, which should be purged with an inert gas. BC-11 (hydrobromide) is soluble in organic solvents such as ethanol and DMSO. It is also soluble in water. The solubility of BC-11 (hydrobromide) in ethanol, DMSO, and water is approximately 20, 25, and 10 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

BC-11 is an inhibitor of the serine protease urokinase-type plasminogen activator (uPA).¹ It reduces cell viability of MDA-MB-231 cells ($EC_{50} = 117 \mu M$).

Reference

1. Longo, A., Librizzi, M., Chuckowree, I.S., et al. Cytotoxicity of the urokinase-plasminogen activator inhibitor carbamimidothioic acid (4-boronophenyl) methyl ester hydrobromide (BC-11) on triple-negative MDA-MB231 breast cancer cells. *Molecules* **20**(6), 9879-9889 (2015).

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, 10/04/2023

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