

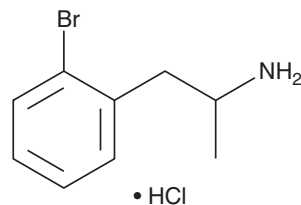
PRODUCT INFORMATION



2-Bromoamphetamine (hydrochloride)

Item No. 9001848

CAS Registry No.: 861006-36-2
Formal Name: 2-bromo- α -methyl-benzeneethanamine, monohydrochloride
Synonyms: 2-BA, o-BA
MF: $C_9H_{12}BrN \cdot HCl$
FW: 250.6
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

4-Bromoamphetamine (4-BA) (Item No. 9001850) is a *para*-substituted amphetamine that acts as a monoamine releasing agent.^{1,2} It is highly neurotoxic, producing long-term depletion of serotonin.^{1,2} 2-BA is a structural isomer of 4-BA, having the bromine at the two rather than the four position. This product is intended only for forensic and research purposes.

References

1. Fuller, R.W., Baker, J.C., Perry, K.W., *et al.* Comparison of 4-chloro-, 4-bromo- and 4-fluoroamphetamine in rats: Drug levels in brain and effects on brain serotonin metabolism. *Neuropharmacology* **14**(10), 739-746 (1975).
2. Harvey, J.A., McMaster, S.E., and Fuller, R.W. Comparison between the neurotoxic and serotonin-depleting effects of various halogenated derivatives of amphetamine in the rat. *J. Pharmacol. Exp. Ther.* **202**(3), 581-589 (1977).

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, 07/12/2023

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