

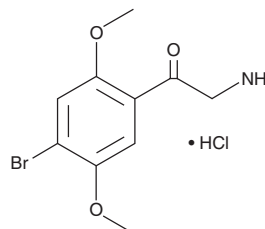
PRODUCT INFORMATION



bk-2C-B (hydrochloride)

Item No. 15328

CAS Registry No.: 2303508-66-7
Formal Name: 2-amino-1-(4-bromo-2,5-dimethoxyphenyl)-ethanone, monohydrochloride
Synonym: bk-4-Bromo-2,5-dimethoxyphenethylamine
MF: C₁₀H₁₂BrNO₃ • HCl
FW: 310.6
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 264, 344 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

A series of 2,5-dimethoxyphenethylamines, substituted at the four position of the aromatic ring, have psychoactive effects.^{1,2} 2C-B (hydrochloride) (Item No. 11734) is a hallucinogenic designer drug that activates the serotonin receptor 5-HT_{2C} (pEC₅₀ = 6.8 for arachidonic acid release) and has emerged as a drug of abuse.¹⁻³ bk-2C-B is a β-ketone analog of 2C-B.⁴ This product is intended for forensic and research purposes.

References

1. Bruno, R., Matthews, A.J., Dunn, M., *et al.* Emerging psychoactive substance use among regular ecstasy users in Australia. *Drug Alcohol Depend.* **124(1-2)**, 19-25 (2012).
2. Moya, P.R., Berg, K.A., Gutiérrez-Hernandez, M.A., *et al.* Functional selectivity of hallucinogenic phenethylamine and phenylisopropylamine derivatives at human 5-hydroxytryptamine (5-HT)_{2A} and 5-HT_{2C} receptors. *J. Pharmacol. Exp. Ther.* **321(3)**, 1054-1061 (2007).
3. Maurer, H.H. Chemistry, pharmacology, and metabolism of emerging drugs of abuse. *Ther. Drug Monit.* **32(5)**, 544-549 (2010).
4. Westphal, F., Junge, T., Girreser, U., *et al.* Structure elucidation of a new designer benzylpiperazine: 4-Bromo-2,5-dimethoxybenzylpiperazine. *Forensic Sci. Int.* **187(1-3)**, 87-96 (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, 06/27/2023

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