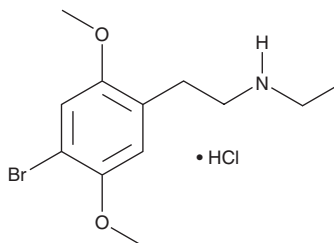


# PRODUCT INFORMATION

## N-ethyl 2C-B (hydrochloride)

Item No. 29387

**CAS Registry No.:** 155638-84-9  
**Formal Name:** 4-bromo-N-ethyl-2,5-dimethoxy-benzeneethanamine, monohydrochloride  
**Synonym:** Ethyl 2C-B  
**MF:** C<sub>12</sub>H<sub>18</sub>BrNO<sub>2</sub> • HCl  
**FW:** 324.6  
**Purity:** ≥98%  
**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

N-ethyl 2C-B (hydrochloride) (Item No. 29387) is an analytical reference standard categorized as a phenethylamine. N-ethyl 2C-B has hallucinogenic properties.<sup>1,2</sup> This product is intended for research and forensic applications.

### References

- Schifano, F., Orsolini, L., Papanti, D., *et al.* NPS: Medical consequences associated with their intake. *Curr. Top. Behav. Neurosci.* **32**, 351-380 (2017).
- Soussan, C., Sundström, K., Andersson, M., *et al.* Poisoning casualties: Alcohol, pharmaceuticals or “legal highs”? Poisoning cases at emergency rooms in the Swedish county Värmland in 2007-2013. *J. Community Med. Health Educ.* **5(6):386** (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, 07/25/2023

#### 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