

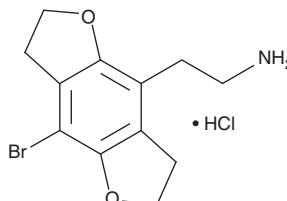
# PRODUCT INFORMATION



## 2C-B-FLY (hydrochloride)

Item No. 11473

**CAS Registry No.:** 178557-21-6  
**Formal Name:** 8-bromo-2,3,6,7-tetrahydrobenzo[1,2-b:4,5-b']difuran-4-ethanamine, monohydrochloride  
**MF:** C<sub>12</sub>H<sub>14</sub>BrNO<sub>2</sub> • HCl  
**FW:** 320.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 203, 315 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Description

A group of phenethylamine derivatives referred to as the FLY compounds, named for their insect-like appearance of two "wing-like" furan or dihydrofuran rings fused on the opposite sides of the central benzene ring, have been identified with allegedly potent hallucinogenic effects.<sup>1</sup> 2C-B-FLY is the dihydrodifuran analog of the Schedule I hallucinogen 4-bromo-2,5-dimethoxyphenethylamine (2C-B).<sup>1,2</sup> It is expected to show similar activity to 2C-B, which acts as a partial agonist at the 5-HT<sub>2A</sub> serotonin receptor and demonstrates high binding affinity for the 5-HT<sub>2B</sub> and 5-HT<sub>2C</sub> serotonin receptors.<sup>3,4</sup> This product is intended for forensic and research purposes.

### References

1. Reed, E.C. and Kiddon, G.S. The characterization of three fly compounds (2C-B-FLY, 3C-B-FLY, and Bromo-Dragon FLY). *DEA Microgram Journal* **5(1-4)**, 1-8 (2007).
2. Pichini, S., Pujadas, M., Marchei, E., et al. Liquid chromatography-atmospheric pressure ionization electrospray mass spectrometry determination of "hallucinogenic designer drugs" in urine of consumers. *J. Pharm. Biomed. Anal.* **47(2)**, 335-342 (2008).
3. Monte, A.P., Marona-Lewicka, D., Parker, M.A., et al. Dihydrobenzofuran analogues of hallucinogens. 3. Models of 4-substituted (2,5-dimethoxyphenyl)alkylamine derivatives with rigidified methoxy groups. *J. Med. Chem.* **39(15)**, 2953-2961 (1996).
4. Chambers, J.J., Kurrasch-Orbaugh, D.M., Parker, M.A., et al. Enantiospecific synthesis and pharmacological evaluation of a series of super-potent, conformationally restricted 5-HT<sub>2A/2C</sub> receptor agonists. *J. Med. Chem.* **44(6)**, 1003-1010 (2001).

#### 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, 12/08/2021

#### 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