

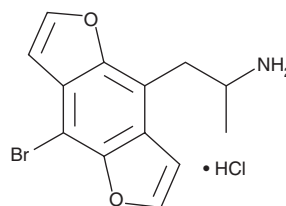
PRODUCT INFORMATION



Bromo-DragonFLY (hydrochloride)

Item No. 11561

CAS Registry No.: 219986-78-4
Formal Name: 8-bromo- α -methyl-benzo[1,2-b:4,5-b']difuran-4-ethanamine, monohydrochloride
Synonyms: 2-Amino-1-(8-bromobenzo[1,2-b:5,4-b']difuran-4-yl)propane, 1-(8-Bromobenzo[1,2-b; 4,5-b']difuran-4-yl)-2-aminopropane, DOB-DFLY
MF: $C_{13}H_{12}BrNO_2 \cdot HCl$
FW: 330.6
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 225, 283, 295 nm
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

Bromo-DragonFLY (hydrochloride) (Item No. 11561) is an analytical reference standard categorized as a phenethylamine.¹ Bromo-DragonFLY has been associated with fatalities and acute overdoses.² It induces the head-twitch response in mice, indicating hallucinogenic potential.³ This product is intended for research and forensic applications.

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). *Microgram J.* **5(1-4)**, 1-8 (2007).
2. Corazza, O., Schifano, F., Farre, M., *et al.* Designer drugs on the internet: A phenomenon out-of-control? The emergence of hallucinogenic drug Bromo-Dragonfly. *Curr. Clin. Pharmacol.* **6(2)**, 125-129 (2011).
3. Halberstadt, A.L., Chatha, M., Stratford, A., *et al.* Comparison of the behavioral responses induced by phenylalkylamine hallucinogens and their tetrahydrobenzodifuran ("FLY") and benzodifuran ("DragonFLY") analogs. *Neuropharmacology* **144**, 368-376 (2019).

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, 07/18/2023

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