

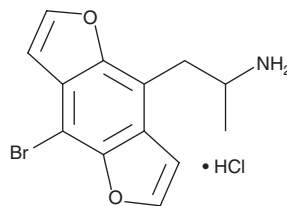
PRODUCT INFORMATION



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
MF: C₁₃H₁₂BrNO₂ • HCl
FW: 330.6
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max} : 225, 283, 295 nm



Laboratory Procedures

For long term storage, we suggest that bromo-dragonFLY (hydrochloride) be stored as supplied at -20°C. It should be stable for at least two years.

Bromo-DragonFLY (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the bromo-dragonFLY (hydrochloride) in the solvent of choice. Bromo-DragonFLY (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of bromo-dragonFLY (hydrochloride) in these solvents is approximately 5 and 1 mg/ml, respectively.

Bromo-DragonFLY (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, bromo-dragonFLY (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Bromo-DragonFLY (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Bromo-DragonFLY is a difuran analog of the Schedule I hallucinogen 4-bromo-2,5-dimethoxyphenethylamine.¹ This designer drug strongly binds serotonin receptors ($K_i = 0.04, 0.19, \text{ and } 0.02 \text{ nM}$ for 5-HT_{2A}, 5-HT_{2B}, and 5-HT_{2C}, respectively).² It is more potent than LSD in drug discrimination studies in rats.² Related to its high potency, bromo-dragonFLY has been associated with fatalities as well as acute intoxications.³ GC-MS, GC-IRD, and solid phase FTIR spectra have been published.¹ This product is intended for research and forensic 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. Parker, M.A., Marona-Lewicka, D., Lucaites, V.L., et al. A novel (benzodifuranyl)aminoalkane with extremely potent activity at the 5-HT_{2A} receptor. *J. Med. Chem.* **41**, 5148-5149 (1998).
3. 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**, 125-129 (2011).

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, 08/27/2015

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