

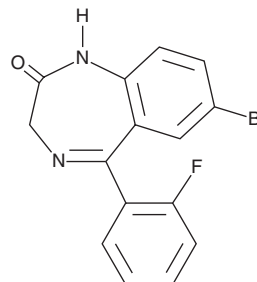
# PRODUCT INFORMATION



## Flubromazepam

Item No. 15157

**CAS Registry No.:** 2647-50-9  
**Formal Name:** 7-bromo-5-(2-fluorophenyl)-1,3-dihydro-2H-1,4-benzodiazepin-2-one  
**MF:** C<sub>15</sub>H<sub>10</sub>BrFN<sub>2</sub>O  
**FW:** 333.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 230, 322 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

Flubromazepam (Item No. 15157) is an analytical reference standard categorized as a benzodiazepine.<sup>1</sup> Flubromazepam has been detected in seized samples and has anxiolytic and sedative properties.<sup>2,3</sup> This product is intended for research and forensic applications.

### References

1. Moosmann, B., Huppertz, L.M., Hutter, M., *et al.* Detection and identification of the designer benzodiazepine flubromazepam and preliminary data on its metabolism and pharmacokinetics. *J. Mass. Spectrom.* **48(11)**, 1150-1159 (2013).
2. Jones, L.E., Stewart, A., Peters, K.L., *et al.* Infrared and raman screening of seized novel psychoactive substances: A large scale study of >200 samples. *Analyst* **141(3)**, 902-909 (2016).
3. Abouchedid, R., Gilks, T., Dargan, P.I., *et al.* Assessment of the availability, cost, and motivations for use over time of the new psychoactive substances-benzodiazepines diclazepam, flubromazepam, and pyrazolam-in the UK. *J. Med. Toxicol.* **14(2)**, 134-143 (2018).

#### 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, 06/26/2023

#### 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