

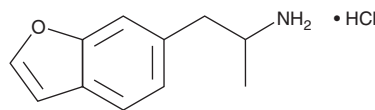
# PRODUCT INFORMATION



## 6-APB (hydrochloride)

Item No. 11079

**CAS Registry No.:** 286834-84-2  
**Formal Name:**  $\alpha$ -methyl-6-benzofuranethanamine, monohydrochloride  
**Synonyms:** 6-(2-Aminopropyl)Benzofuran, Benzo Fury  
**MF:**  $C_{11}H_{13}NO \cdot HCl$   
**FW:** 211.7  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 207, 246, 276, 285 nm  
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 2$  years



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

### Description

6-APB (hydrochloride) (Item No. 11079) is an analytical reference standard categorized as a benzofuran. 6-APB is a metabolite of 6-MAPB (Item No. 14110).<sup>1</sup> It is a potential metabolite of 6-EAPB (Item No. 9002053). 6-APB has been found in seized samples suspected of containing novel psychoactive substances (NPS).<sup>2</sup> This product is intended for research and forensic applications.

### References

1. Welter, J., Brandt, S.D., Kavanagh, P., *et al.* Metabolic fate, mass spectral fragmentation, detectability, and differentiation in urine of the benzofuran designer drugs 6-APB and 6-MAPB in comparison to their 5-isomers using GC-MS and LC-(HR)-MS<sup>n</sup> techniques. *Anal. Bioanal. Chem.* **407(12)**, 3457-3470 (2012).
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).

#### 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, 12/09/2022

#### 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