# **PRODUCT INFORMATION**



# 4-Bromomethcathinone (hydrochloride)

Item No. 12089

CAS Registry No.: 135333-27-6

Formal Name: 1-(4-bromophenyl)-2-

(methylamino)-1-propanone,

monohydrochloride

Synonyms: 4-BMC, Brephedrone MF: C<sub>10</sub>H<sub>12</sub>BrNO • HCl

FW: 278.6 **Purity:** ≥98% Supplied as: A neat solid Storage: -20°C Stability: ≥2 years

HCl

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

## Description

4-Bromomethcathinone (hydrochloride) (Item No. 12089) is an analytical reference material categorized as a cathinone. 4-Bromomethcathinone has been detected in capsules seized by law enforcement. 1 This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

### Reference

1. Machado, Y., Neto, J.C., Barbosa, P.E.N., et al. Brephedrone: A new psychoactive substance seized in Brazil. Forensic Sci. Int. 275, 302-307 (2017).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information Buyer agrees to purchase the material can be found on our website.

Copyright Cayman Chemical Company, 08/31/2020

# **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