

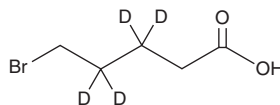
PRODUCT INFORMATION



5-Bromopentanoic Acid-d₄

Item No. 30913

CAS Registry No.: 42932-62-7
Formal Name: 5-bromo-pentanoic-3,3,4,4-d₄ acid
Synonyms: 5-Bromovaleric Acid-d₄, δ-Bromovaleric Acid-d₄
MF: C₅H₅BrD₄O₂
FW: 185.1
Chemical Purity: ≥98% (5-Bromopentanoic Acid)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

5-Bromopentanoic acid-d₄ is intended for use as an internal standard for the quantification of 5-bromopentanoic acid by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

5-Bromopentanoic acid-d₄ is supplied as a solid. A stock solution may be made by dissolving the 5-bromopentanoic acid-d₄ in the solvent of choice, which should be purged with an inert gas. 5-Bromopentanoic acid-d₄ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 5-bromopentanoic acid-d₄ in these solvents is approximately 30 mg/ml.

Description

5-Bromopentanoic acid is a synthetic intermediate.¹ It has been used in the synthesis of PI3K inhibitors.

Reference

1. Li, Y., Zeng, Q., Wang, R., *et al.* Synthesis and discovery of mitochondria-targeting oleanolic acid derivatives for potential PI3K inhibition. *Fitoterapia* **162**, 105291 (2022).

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, 04/05/2023

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