

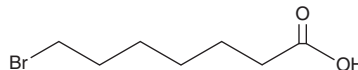
PRODUCT INFORMATION



7-Bromoheptanoic Acid

Item No. 30420

CAS Registry No.: 30515-28-7
MF: $C_7H_{13}BrO_2$
FW: 209.1
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

7-Bromoheptanoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 7-bromoheptanoic acid in the solvent of choice, which should be purged with an inert gas. 7-Bromoheptanoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 7-bromoheptanoic acid in these solvents is approximately 30 mg/ml.

7-Bromoheptanoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 7-bromoheptanoic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 7-Bromoheptanoic acid has a solubility of approximately 0.16 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

7-Bromoheptanoic acid is a building block.^{1,2} It has been used in the synthesis of azide-based nicotinamide phosphoribosyltransferase (Nampt) inhibitors with anticancer activity and SAHA (Item No. 10009929) derivatives that inhibit histone deacetylases (HDACs).

References

1. Colombano, G., Travelli, C., Galli, U., *et al.* A novel potent nicotinamide phosphoribosyltransferase inhibitor synthesized via click chemistry. *J. Med. Chem.* **53**(2), 616-623 (2010).
2. Suzuki, T., Nagano, Y., Kouketsu, A., *et al.* Novel inhibitors of human histone deacetylases: Design, synthesis, enzyme inhibition, and cancer cell growth inhibition of SAHA-based non-hydroxamates. *J. Med. Chem.* **48**(4), 1019-1032 (2005).

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, 11/18/2022

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