

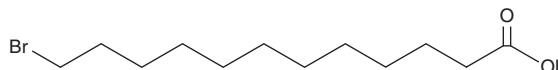
# PRODUCT INFORMATION



## 12-Bromododecanoic Acid

Item No. 30465

CAS Registry No.: 73367-80-3  
Synonyms: 12-Bromo-C12:0,  
12-Bromododecanoate,  
12-Bromolauric Acid, NSC 660375  
MF:  $C_{12}H_{23}BrO_2$   
FW: 279.2  
Purity:  $\geq 95\%$   
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

12-Bromododecanoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 12-bromododecanoic acid in the solvent of choice, which should be purged with an inert gas. 12-Bromododecanoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 12-bromododecanoic acid in these solvents is approximately 30 mg/ml.

### Description

12-Bromododecanoic acid is a halogenated form of lauric acid (Item No. 10006626). It has been used in the synthesis of clickable forms of myristic acid (Item No. 13351) and as a model fatty acid ligand to determine the X-ray crystal structure of bovine  $\beta$ -lactoglobulin-ligand complexes.<sup>1,2</sup> 12-Bromododecanoic acid (10  $\mu$ g/ml) reduces virion DNA in the culture supernatant of primary hepatocytes isolated from a duckling model of hepatitis B virus (HBV) infection and inhibits HIV replication in CEM-SS T cells ( $EC_{50} = 38 \mu M$ ).<sup>3</sup>

### References

1. Heal, W.P., Wickramasinghe, S.R., Leatherbarrow, R.J., et al. N-Myristoyl transferase-mediated protein labelling *in vivo*. *Org. Biomol. Chem.* **6**(13), 2308-2315 (2008).
2. Qin, B.Y., Creamer, L.K., Baker, E.N., et al. 12-Bromododecanoic acid binds inside the calyx of bovine  $\beta$ -lactoglobulin. *FEBS Lett.* **438**(3), 272-278 (1998).
3. Parang, K., Wiebe, L.I., Knaus, E.E., et al. In vitro antiviral activities of myristic acid analogs against human immunodeficiency and hepatitis B viruses. *Antiviral Res.* **34**(3), 75-90 (1997).

#### 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, 12/15/2022

#### 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