

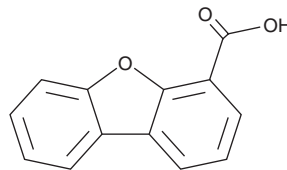
PRODUCT INFORMATION



Dibenzofuran-4-carboxylic Acid

Item No. 36681

CAS Registry No.: 2786-05-2
Formal Name: 4-dibenzofurancarboxylic acid
Synonym: NSC 190611
MF: C₁₃H₈O₃
FW: 212.2
Purity: ≥95%
UV/Vis.: λ_{max}: 217, 241, 291 nm
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

Dibenzofuran-4-carboxylic acid is supplied as a solid. A stock solution may be made by dissolving the dibenzofuran-4-carboxylic acid in the solvent of choice, which should be purged with an inert gas. Dibenzofuran-4-carboxylic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of dibenzofuran-4-carboxylic acid in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

Dibenzofuran-4-carboxylic acid is a synthetic precursor.^{1,2} It has been used in the synthesis of light-emitting polycyclic aromatic hydrocarbons (PAHs) and *S. mansoni* histone deacetylase 8 (HDAC8) inhibitors.

References

1. Molotkov, A.P., Arsenov, M.A., Kapustin, D.A., *et al.* Effect of Cp-ligand methylation on rhodium(III)-catalyzed annulations of aromatic carboxylic acids with alkynes: Synthesis of isocoumarins and PAHs for organic light-emitting devices. *Chempluschem* **85(2)**, 334-345 (2020).
2. Ghazy, E., Heimburg, T., Lancelot, J., *et al.* Synthesis, structure-activity relationships, cocrystallization and cellular characterization of novel smHDAC8 inhibitors for the treatment of schistosomiasis. *Eur. J. Med. Chem.* **225:113745**, (2021).

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, 06/16/2023

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