

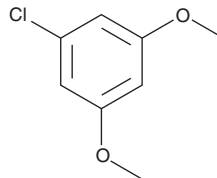
PRODUCT INFORMATION



5-Chlororesorcinol dimethyl ether

Item No. 10007127

CAS Registry No.: 7051-16-3
Formal Name: 1-chloro-3,5-dimethoxy-benzene
Synonym: NSC 220095
MF: C₈H₉ClO₂
FW: 172.6
Purity: ≥98%
Supplied as: A liquid
Storage: Room Temperature
Stability: ≥2 years



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

Laboratory Procedures

5-Chlororesorcinol dimethyl ether is supplied as a liquid. A stock solution may be made by dissolving the 5-chlororesorcinol dimethyl ether in the solvent of choice. 5-Chlororesorcinol dimethyl ether is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 5-chlororesorcinol dimethyl ether in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

5-chlororesorcinol dimethyl ether is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the solution of 5-chlororesorcinol dimethyl ether should be diluted with the aqueous buffer of choice. 5-Chlororesorcinol dimethyl ether has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method.

Description

5-Chlororesorcinol dimethyl ether is a synthetic intermediate useful for pharmaceutical synthesis.

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, 06/07/2018

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