

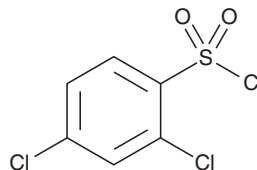
# PRODUCT INFORMATION



## 2,4-Dichlorobenzenesulfonyl Chloride

Item No. 30450

CAS Registry No.: 16271-33-3  
Formal Name: 2,4-dichloro-benzenesulfonyl chloride  
MF:  $C_6H_3Cl_3O_2S$   
FW: 245.5  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{max}$ : 246 nm  
Supplied as: A 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

2,4-Dichlorobenzenesulfonyl chloride is supplied as a solid. A stock solution may be made by dissolving the 2,4-dichlorobenzenesulfonyl chloride in the solvent of choice, which should be purged with an inert gas. 2,4-Dichlorobenzenesulfonyl chloride is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2,4-dichlorobenzenesulfonyl chloride in these solvents is approximately 25, 10, and 15 mg/ml, respectively.

2,4-Dichlorobenzenesulfonyl chloride is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2,4-dichlorobenzenesulfonyl chloride should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 2,4-Dichlorobenzenesulfonyl chloride has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

2,4-Dichlorobenzenesulfonyl chloride is a heterocyclic building block.<sup>1</sup> It has been used in the synthesis of nicotinyl arylsulfonylhydrazides that have anticancer and antibacterial activities *in vitro*.

### Reference

1. Kamal, A., Khan, M.N.A., Reddy, K.S., *et al.* Synthesis of a new class of 2-anilino substituted nicotinyl arylsulfonylhydrazides as potential anticancer and antibacterial agents. *Bioorg. Med. Chem.* **15**(2), 1004-1013 (2007).

#### 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, 11/02/2022

#### 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