

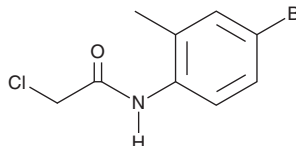
PRODUCT INFORMATION



Cosan-528

Item No. 37537

CAS Registry No.: 96686-51-0
Formal Name: N-(4-bromo-2-methylphenyl)-2-chloro-acetamide
MF: $C_9H_9BrClNO$
FW: 262.5
Purity: $\geq 98\%$
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

Cosan-528 is supplied as a solid. A stock solution may be made by dissolving the cosan-528 in the solvent of choice, which should be purged with an inert gas. Cosan-528 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of cosan-528 in DMF is approximately 25 mg/ml and approximately 10 mg/ml in ethanol and DMSO.

Description

Cosan-528 is a fungicide.¹ It is active against the fungus *A. pullulans* (MIC = 8 ppm). Cosan-528 completely inhibits the growth of *A. pullulans*, *A. niger*, and *Penicillium* on paint-coated white pine sapwood for at least six weeks.

Reference

1. Gagliani, K. and Eilender, A.L. Industrial fungicides. *Cosan Chemical Corporation*. 0137729A2 (1985).

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, 06/06/2023

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