

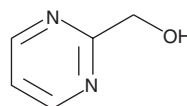
# PRODUCT INFORMATION



## 2-(hydroxymethyl)-Pyrimidine

Item No. 10007112

**CAS Registry No.:** 42839-09-8  
**Formal Name:** 2-pyrimidinemethanol  
**MF:** C<sub>5</sub>H<sub>6</sub>N<sub>2</sub>O  
**FW:** 110.1  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 248, 286 nm nm  
**Supplied as:** A crystalline 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

2-(Hydroxymethyl)-pyrimidine is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-(hydroxymethyl)-pyrimidine in the solvent of choice, which should be purged with an inert gas. 2-(Hydroxymethyl)-Pyrimidine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2-(hydroxymethyl)-pyrimidine in ethanol is approximately 100 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

2-(hydroxymethyl)-Pyrimidine 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, 12/11/2022

#### 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