

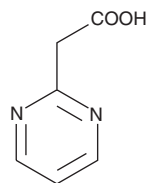
# PRODUCT INFORMATION



## 2-Pyrimidine Acetic Acid

Item No. 10007116

CAS Registry No.: 66621-73-6  
Formal Name: 2-pyrimidineacetic acid  
MF:  $C_6H_6N_2O_2$   
FW: 138.1  
Purity:  $\geq 95\%$   
UV/Vis.:  $\lambda_{max}$ : 250 nm  
Supplied as: A crystalline 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-Pyrimidine acetic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-pyrimidine acetic acid in the organic solvent of choice, which should be purged with an inert gas. 2-Pyrimidine acetic acid is soluble in the organic solvent dimethyl formamide at a concentration of approximately 1 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 2-pyrimidine acetic acid can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of 2-pyrimidine acetic acid in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

2-Pyrimidine acetic acid 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