

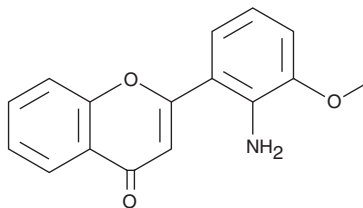
PRODUCT INFORMATION



2-methyl-1,3-Cyclohexanedione

Catalog No. 10007138

CAS Registry No.: 1193-55-1
Formal Name: 2-methyl-1,3-cyclohexanedione
Synonym: NSC 54460
MF: C₇H₁₀O₂
FW: 126.2
Purity: ≥95%
UV/Vis.: λ_{max}: 262 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-methyl-1,3-Cyclohexanedione is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-methyl-1,3-cyclohexanedione in the solvent of choice, which should be purged with an inert gas. 2-methyl-1,3-Cyclohexanedione is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2-methyl-1,3-cyclohexanedione in ethanol is approximately 2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

2-methyl-1,3-Cyclohexanedione is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2-methyl-1,3-cyclohexanedione should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2-methyl-1,3-Cyclohexanedione has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

2-methyl-1,3-Cyclohexanedione 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, 12/12/2022

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