

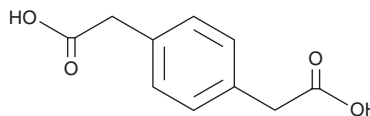
PRODUCT INFORMATION



1,4-Phenylenediacetic Acid

Item No. 30449

CAS Registry No.: 7325-46-4
Formal Name: 1,4-benzenediacetic acid
Synonyms: *p*-Benzenediacetic Acid,
p-Phenylenediacetic Acid
MF: C₁₀H₁₀O₄
FW: 194.2
Purity: ≥98%
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

1,4-Phenylenediacetic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,4-phenylenediacetic acid in the solvent of choice, which should be purged with an inert gas. 1,4-Phenylenediacetic acid is soluble in ethanol and dimethyl formamide. The solubility of 1,4-phenylenediacetic acid in these solvents is approximately 30 and 5 mg/ml, respectively.

Description

1,4-Phenylenediacetic acid is a building block.^{1,2} It has been used in the synthesis of metal-ligand coordination polymers.

References

1. Braverman, M.A. and LaDuca, R.L. Luminescent two- and three-dimensional zinc coordination polymers containing isomers of phenylenediacetate and a kinked tethering organodiimine. *Cryst. Growth Des.* **7**(11), 2343-2351 (2007).
2. Yang, G.-P., Wang, Y.-Y., Zhang, W.-H., *et al.* A series of Zn(II) coordination complexes derived from isomeric phenylenediacetic acid and dipyriddy ligands: syntheses, crystal structures, and characterizations. *CrystEngComm* **12**(5), 1509-1517 (2010).

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, 10/27/2022

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