

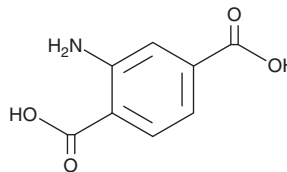
PRODUCT INFORMATION



2-Aminoterephthalic Acid

Item No. 45551

CAS Registry No.: 10312-55-7
Formal Name: 2-amino-1,4-benzenedicarboxylic acid
Synonym: 2-ATA
MF: C₈H₇NO₄
FW: 181.1
Purity: ≥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

2-Aminoterephthalic acid is supplied as a solid. A stock solution may be made by dissolving the 2-aminoterephthalic acid in the solvent of choice, which should be purged with an inert gas. 2-Aminoterephthalic acid is slightly soluble (0.1-1 mg/ml) in methanol.

Description

2-Aminoterephthalic acid is a building block that has been used as a ligand in coordination chemistry and in the synthesis of metal-organic frameworks (MOFs).^{1,2}

References

1. Yang, Y., Lin, R., Ge, L., *et al.* Synthesis and characterization of three amino-functionalized metal-organic frameworks based on the 2-aminoterephthalic ligand. *Dalton Trans.* **44**(17), 8190-8197 (2015).
2. Arias-Pardilla, J., Salavagione, H.J., Barbero, C., *et al.* Study of the chemical copolymerization of 2-aminoterephthalic acid and aniline.: Synthesis and copolymer properties. *Eur. Polym. J.* **42**(7), 1521-1532 (2006).

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, 04/30/2026

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