

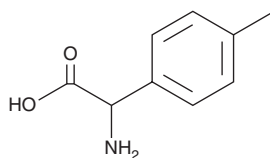
PRODUCT INFORMATION



2-Amino-2-(*p*-tolyl)acetic Acid

Item No. 42294

CAS Registry No.: 13227-01-5
Formal Name: α -amino-4-methyl-benzeneacetic acid
Synonyms: 2-Amino-2-(*para*-tolyl)acetic Acid,
2-Amino-2-(4-tolyl)acetic Acid,
4-Methylphenyl Glycine
MF: $C_9H_{11}NO_2$
FW: 165.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

2-Amino-2-(*p*-tolyl)acetic acid is supplied as a solid. A stock solution may be made by dissolving the 2-amino-2-(*p*-tolyl)acetic acid in the solvent of choice, which should be purged with an inert gas. 2-Amino-2-(*p*-tolyl)acetic acid is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

2-Amino-2-(*p*-tolyl)acetic acid is a synthetic intermediate.¹ It has been used in the synthesis of phenylalanine ammonia-lyase inhibitors.

Reference

1. Zoń, J., Amfrhein, N., and Gancarz, R. Inhibitors of phenylalanine ammonia-lyase: 1-Aminobenzylphosphonic acids substituted in the benzene ring. *Phytochemistry* **59**(1), 9-21 (2002).

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, 12/11/2024

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