

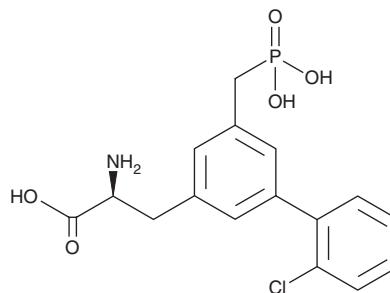
PRODUCT INFORMATION



SDZ 220-581

Item No. 33344

CAS Registry No.: 174575-17-8
Formal Name: αS-amino-2'-chloro-5-(phosphonomethyl)-[1,1'-biphenyl]-3-propanoic acid
MF: C₁₆H₁₇ClNO₅P
FW: 369.7
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

SDZ 220-581 is supplied as a solid. A stock solution may be made by dissolving the SDZ 220-581 in the solvent of choice, which should be purged with an inert gas. SDZ 220-581 is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

Description

SDZ 220-581 is an NMDA receptor antagonist (apparent $pA_2 = 7.8$).¹ It reduces the number of convulsions induced by maximal electroshock (MES) in mice ($ED_{50} = <3.2$ mg/kg).² SDZ 220-581 (1.25 mg/kg) decreases lesion area in a mouse model of cerebral ischemia induced by middle cerebral artery occlusion (MCAO). It reduces mechanical hyperalgesia induced by complete Freund's adjuvant in mice when administered at doses of 10 or 30 mg/kg.

References

1. Müller, W., Kipfer, P., Lowe, D.A., *et al.* Syntheses of biphenyl analogues of AP7, a new class of competitive N-methyl-D-aspartate (NMDA) receptor antagonists. *Helv. Chim. Acta* **78**, 2026-2035 (1995).
2. Urwyler, S., Campbell, E., Fricker, G., *et al.* Biphenyl-derivatives of 2-amino-7-phosphono-heptanoic acid, a novel class of potent competitive N-methyl-D-aspartate receptor antagonists--II. Pharmacological characterization *in vivo*. *Neuropharmacology* **35(6)**, 655-669 (1996).

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, 08/11/2025

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