

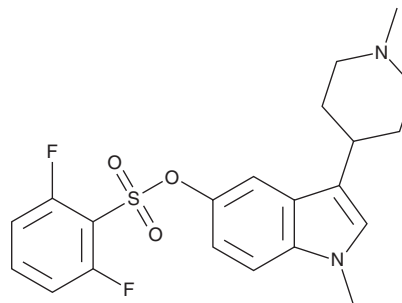
# PRODUCT INFORMATION



**SDZ 220-040**

Item No. 43672

**CAS Registry No.:** 174575-40-7  
**Formal Name:** αS-amino-2',4'-dichloro-4-hydroxy-5-(phosphonomethyl)-[1,1'-biphenyl]-3-propanoic acid  
**MF:** C<sub>16</sub>H<sub>16</sub>Cl<sub>2</sub>NO<sub>6</sub>P  
**FW:** 420.2  
**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 220-040 is supplied as a solid. A stock solution may be made by dissolving the SDZ 220-040 in the solvent of choice, which should be purged with an inert gas. SDZ 220-040 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

SDZ 220-040 is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

## Description

SDZ 220-040 is an NMDA receptor antagonist (apparent  $pA_2 = 8.75$ ).<sup>1</sup> It reduces the number of convulsions induced by maximal electroshock (MES) in mice ( $ED_{50} = 3.2-10$  mg/kg).<sup>2</sup>

## 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, 07/17/2025

## 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](http://WWW.CAYMANCHEM.COM)