

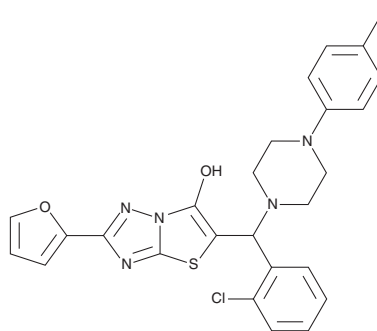
# PRODUCT INFORMATION



## GluN1/3A-IN-2

Item No. 45520

**CAS Registry No.:** 887220-24-8  
**Formal Name:** 5-[(2-chlorophenyl)[4-(4-fluorophenyl)-1-piperazinyl]methyl]-2-(2-furanyl)-thiazolo[3,2-b][1,2,4]triazol-6-ol  
**Synonyms:** Compound DD-1, , GluN1/3A Inhibitor 2  
**MF:** C<sub>25</sub>H<sub>21</sub>ClFN<sub>5</sub>O<sub>2</sub>S  
**FW:** 510.0  
**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

GluN1/3A-IN-2 is supplied as a solid. A stock solution may be made by dissolving the GluN1/3A-IN-2 in the solvent of choice, which should be purged with an inert gas. GluN1/3A-IN-2 is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

GluN1/3A-IN-2 is an inhibitor of NR1/3A and NR1/3B subunit-containing NMDA receptors ( $IC_{50}$ s = 1.56 and 0.12  $\mu$ M, respectively).<sup>1</sup> It is selective for NR1/3A and NR1/3B subunit-containing NMDA receptors over those containing NR1/2A subunits ( $IC_{50}$  = >30  $\mu$ M). GluN1/3A-IN-2 inhibits glycine-induced currents in a whole-cell patch-clamp assay using HEK293 cells expressing NR1/3A subunit-containing NMDA receptors.

### Reference

1. Li, S.W., Zeng, Y., Wu, S.N., *et al.* Discovery of selective GluN1/GluN3A NMDA receptor inhibitors using integrated AI and physics-based approaches. *Acta. Pharmacol. Sin.* **47(1)**, 22-31 (2026).

#### 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, 04/24/2026

#### 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