

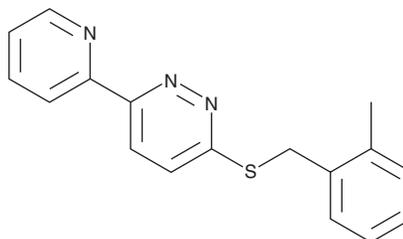
# PRODUCT INFORMATION



**LDN-0212320**

Item No. 21650

**CAS Registry No.:** 894002-50-7  
**Formal Name:** 3-[[[(2-methylphenyl)methyl]thio]-6-(2-pyridinyl)-pyridazine  
**Synonym:** OSU-0212320  
**MF:** C<sub>17</sub>H<sub>15</sub>N<sub>3</sub>S  
**FW:** 293.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 294 nm  
**Supplied as:** A crystalline 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

LDN-0212320 is supplied as a crystalline solid. A stock solution may be made by dissolving the LDN-0212320 in the solvent of choice, which should be purged with an inert gas. LDN-0212320 is soluble in organic solvents such as ethanol and DMSO. The solubility of LDN-0212320 in these solvents is approximately 5 and 25 mg/ml, respectively.

## Description

LDN-0212320 is an activator of excitatory amino acid transporter 2 (EAAT2) that increases expression of EAAT2 in PA-EAAT2 cells in a concentration- and time-dependent manner (EC<sub>50</sub> = 1.83 μM).<sup>1</sup> It selectively increases EAAT2 protein levels over EAAT1 and EAAT3 protein levels and increases glutamate uptake in primary dissociated neuron and astrocyte mixed cultures when used at concentrations ranging from 1 to 3 μM. LDN-0212320 (3 μM) decreases neuronal loss and degeneration in primary dissociated neuron and astrocyte mixed cultures. *In vivo*, LDN-0212320 (40 mg/kg, i.p.) increases EAAT2 protein levels and glutamate uptake in mouse forebrain. It also delays motor function decline and increases lifespan in the SOD1G93A transgenic mouse model of amyotrophic lateral sclerosis (ALS).

## Reference

1. Kong, Q., Chang, L.-C., Takahashi, K., *et al.* Small-molecule activator of glutamate transporter EAAT2 translation provides neuroprotection. *J. Clin. Invest.* **124**(3), 1255-1267 (2014).

### 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, 11/14/2022

## 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