

# PRODUCT INFORMATION



**LY393558**

Item No. 32775

**CAS Registry No.:** 271780-64-4  
**Formal Name:** 1-[2-[4-(6-fluoro-1H-indol-3-yl)-3,6-dihydro-1(2H)-pyridinyl]ethyl]-3,4-dihydro-3-(1-methylethyl)-6-(methylsulfonyl)-1H-2,1,3-benzothiadiazine, 2,2-dioxide

**MF:** C<sub>26</sub>H<sub>31</sub>FN<sub>4</sub>O<sub>4</sub>S<sub>2</sub>

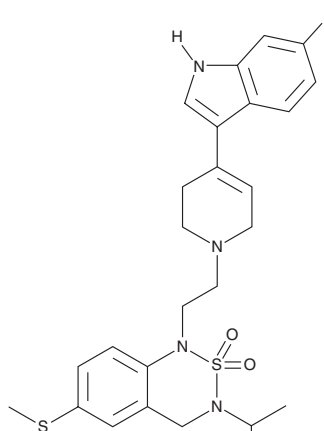
**FW:** 546.7

**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

LY393558 is supplied as a solid. A stock solution may be made by dissolving the LY393558 in the solvent of choice, which should be purged with an inert gas. LY393558 is soluble in organic solvents such as ethanol and DMSO. The solubility of LY393558 in these solvents is approximately 10 mg/ml.

## Description

LY393558 is an inhibitor of serotonin (5-HT) reuptake ( $IC_{50} = 3.3$  nM in rat cortical synaptosomes).<sup>1</sup> It is selective for serotonin reuptake over norepinephrine reuptake in rat cortical synaptosomes ( $IC_{50} = 2,510$  nM) but does inhibit 5-HT-induced [<sup>35</sup>S]GTPγS binding in cells expressing human 5-HT<sub>1B</sub> or 5-HT<sub>1D</sub> receptors ( $K_{iS} = 0.89$  and 10.4 nM, respectively). LY393558 (30 mg/kg per day) reduces pulmonary vascular remodeling, increases in right ventricular systolic pressure, and the development of right ventricular hypertrophy in a mouse model of pulmonary arterial hypertension induced by overexpression of the serotonin transporter and hypoxia.<sup>2</sup>

## References

1. Pullar, I.A., Boot, J.R., Carney, S.L., *et al.* In vitro activity of LY393558, an inhibitor of the 5-hydroxytryptamine transporter with 5-HT<sub>1B/1D/2</sub> receptor antagonist properties. *Eur. J. Pharmacol.* **432**(1), 9-17 (2001).
2. Morecroft, I., Pang, L., Baranowska, M., *et al.* In vivo effects of a combined 5-HT<sub>1B</sub> receptor/SERT antagonist in experimental pulmonary hypertension. *Cardiovasc. Res.* **85**(3), 593-603 (2010).

### 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, 10/26/2022

## 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