

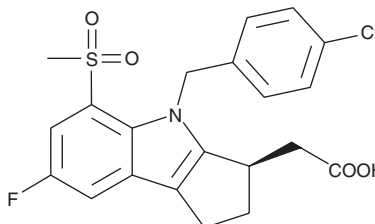
# PRODUCT INFORMATION



**MK-0524**

Item No. 10009835

**CAS Registry No.:** 571170-77-9  
**Formal Name:** (3R)-4-[(4-chlorophenyl)methyl]-7-fluoro-1,2,3,4-tetrahydro-5-(methylsulfonyl)-cyclopent[b]indole-3-acetic acid  
**Synonym:** Laropiprant  
**MF:** C<sub>21</sub>H<sub>19</sub>ClFNO<sub>4</sub>S  
**FW:** 435.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 243, 311 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

MK-0524 is supplied as a crystalline solid. A stock solution may be made by dissolving the MK-0524 in the solvent of choice, which should be purged with an inert gas. MK-0524 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MK-0524 in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

MK-0524 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MK-0524 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MK-0524 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

MK-0524 is a potent, selective prostaglandin D<sub>2</sub> (PGD<sub>2</sub>) receptor DP<sub>1</sub> antagonist with K<sub>i</sub> values of 0.57 nM and 0.75 μM for the DP<sub>1</sub> and DP<sub>2</sub> receptors, respectively.<sup>1</sup> It inhibits PGD<sub>2</sub>-induced accumulation of cAMP in both washed platelets and platelet-rich plasma with IC<sub>50</sub> values of 0.09 and 4.0 nM, respectively.<sup>1</sup> In a sheep model of allergic rhinitis, 0.1 mg/kg MK-0524 completely blocked PGD<sub>2</sub>-induced nasal congestion.<sup>1</sup> At a dose of 4 mg/kg, MK-0524 suppressed a nicotinic acid-induced vasodilatory response by 80% in a mouse model of flushing, an undesirable side-effect of niacin treatment for dyslipidemia.<sup>2</sup> In an *in vitro* model of amyotrophic lateral sclerosis, MK-0524 has been shown to partially protect cultured motor neurons from PGD<sub>2</sub>-induced toxicity.<sup>3,4</sup>

## References

1. Sturino, C.F., O'Neill, G., Lachance, N., *et al.* Discovery of a potent and selective prostaglandin D<sub>2</sub> receptor antagonist, [(3R)-4-(4-chloro-benzyl)-7-fluoro-5-(methylsulfonyl)-1,2,3,4-tetrahydrocyclopent[b]indol-3-yl]-acetic acid. *J. Med. Chem.* **50**(4), 794-806 (2007).
2. Cheng, K., Wu, T.-J., Wu, K.K., *et al.* Antagonism of the prostaglandin D<sub>2</sub> receptor 1 suppresses nicotinic acid-induced vasodilation in mice and humans. *Proc. Natl. Acad. Sci. USA* **103**(17), 6682-6687 (2006).
3. Marchetto, M.C.N., Muotri, A.R., Mu, Y., *et al.* Non-cell-autonomous effect of human SOD1G37R astrocytes on motor neurons derived from human embryonic stem cells. *Cell. Stem Cell* **3**(6), 649-657 (2008).
4. Di Giorgio, F.P., Boulting, G.L., Bobrowicz, S., *et al.* Human embryonic stem cell-derived motor neurons are sensitive to the toxic effect of glial cells carrying an ALS-causing mutation. *Cell. Stem Cell.* **3**(6), 637-648 (2008).

### 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, 12/06/2022

## 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