

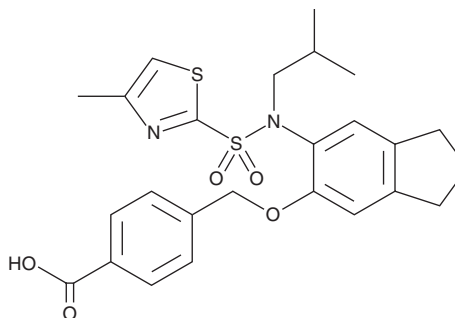
PRODUCT INFORMATION



ONO-8130

Item No. 19118

CAS Registry No.: 459841-96-4
Formal Name: 4-[[[2,3-dihydro-6-[(2-methylpropyl) [(4-methyl-2-thiazolyl)sulfonyl] amino]-1H-inden-5-yl]oxy]methyl]-benzoic acid
MF: C₂₅H₂₈N₂O₅S₂
FW: 500.6
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 280 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

ONO-8130 is supplied as a crystalline solid. A stock solution may be made by dissolving the ONO-8130 in the solvent of choice, which should be purged with an inert gas. ONO-8130 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ONO-8130 in these solvents is approximately 2, 3, and 1 mg/ml, respectively.

ONO-8130 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ONO-8130 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ONO-8130 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

ONO-8130 is an orally bioavailable antagonist of the prostaglandin E₂ (PGE₂) receptor EP₁ (K_i = 1.9 nM)^{1,2} It exerts a more than 1,000-fold selectivity for EP₁ compared with other EP receptors.² ONO-8130 suppresses pain associated with cyclophosphamide-induced cystitis in mice at 0.3-30 mg/kg.¹ It also blocks PGE₂-induced contraction and eliminates spontaneous tone of guinea pig trachea tested in tissue organ baths.²

References

1. Miki, T., Matsunami, M., Nakamura, S., *et al.* ONO-8130, a selective prostanoid EP₁ receptor antagonist, relieves bladder pain in mice with cyclophosphamide-induced cystitis. *Pain* **152**(6), 1373-1381 (2011).
2. S  fholm, J., Dahl  n, S.E., Delin, I., *et al.* PGE₂ maintains the tone of the guinea pig trachea through a balance between activation of contractile EP₁ receptors and relaxant EP₂ receptors. *Br. J. Pharmacol.* **168**, 794-806 (2013).

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, 01/03/2023

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