

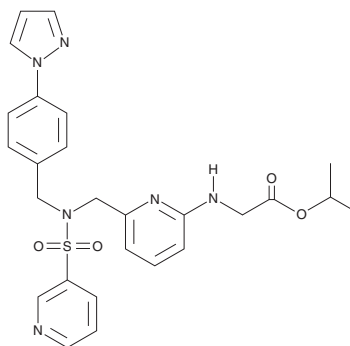
PRODUCT INFORMATION



Omidenepag Isopropyl

Item No. 41046

CAS Registry No.: 1187451-19-9
Formal Name: N-[6-[[[4-(1H-pyrazol-1-yl)phenyl]methyl](3-sulfonyl)amino]methyl]-2-pyridinyl]-glycine, 1-methylethyl ester
Synonyms: DE 11, OMDI
MF: C₂₆H₂₈N₆O₄S
FW: 520.6
Purity: ≥95%
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

Omidenepag isopropyl is supplied as a solid. A stock solution may be made by dissolving the omidenepag isopropyl in the solvent of choice, which should be purged with an inert gas. Omidenepag isopropyl is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Omidenepag isopropyl is a prodrug form of omidenepag (Item No. 41045), a non-prostaglandin agonist of the prostaglandin E₂ (PGE₂) receptor subtype EP₂.¹ It reduces intraocular pressure (IOP) in ocular normotensive and laser-induced ocular hypertensive cynomolgus monkeys when topically administered at a concentration of 0.01%/eye.^{1,2} Formulations containing omidenepag isopropyl have been used in the treatment of open-angle glaucoma and ocular hypertension.

References

1. Iwamura, R., Tanaka, M., Okanari, E., *et al.* Identification of a selective, non-prostanoid EP₂ receptor agonist for the treatment of glaucoma: Omidenepag and its prodrug omidenepag isopropyl. *J. Med. Chem.* **61**(15), 6869-6891 (2018).
2. Kiriara, T., Taniguchi, T., Yamamura, K., *et al.* Pharmacologic characterization of omidenepag isopropyl, a novel selective EP₂ receptor agonist, as an ocular hypotensive agent. *Invest. Ophthalmol. Vis. Sci.* **59**(1), 145-153 (2018).

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, 10/24/2024

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