

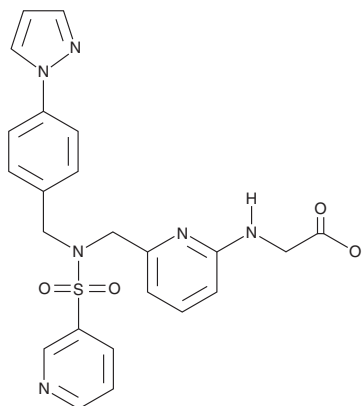
# PRODUCT INFORMATION



## Omidenepag (free acid)

Item No. 41045

**CAS Registry No.:** 1187451-41-7  
**Formal Name:** N-[6-[[[4-(1H-pyrazol-1-yl)phenyl]methyl](3-pyridinylsulfonyl)amino]methyl]-2-pyridinyl]-glycine  
**Synonym:** UR-7276  
**MF:** C<sub>23</sub>H<sub>22</sub>N<sub>6</sub>O<sub>4</sub>S  
**FW:** 478.5  
**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

Omidenepag (free acid) is supplied as a solid. A stock solution may be made by dissolving the omidenepag (free acid) in the solvent of choice, which should be purged with an inert gas. Omidenepag (free acid) is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

Omidenepag is a non-prostaglandin agonist of the prostaglandin E<sub>2</sub> (PGE<sub>2</sub>) receptor subtype EP<sub>2</sub> and an active metabolite of the prodrug omidenepag isopropyl ester.<sup>1</sup> It selectively binds to human EP<sub>2</sub> over EP<sub>1</sub>, EP<sub>3</sub>, and EP<sub>4</sub> receptors (IC<sub>50</sub>s = 10, >10,000, >10,000, and 5,480 nM, respectively) and induces cAMP accumulation in HEK293 cells expressing human EP<sub>2</sub> receptors (EC<sub>50</sub> = 1.1 nM).

### 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).

#### 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, 10/16/2024

#### 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