

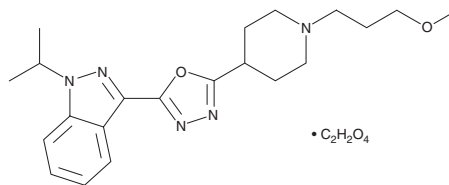
PRODUCT INFORMATION



Usmarapride (oxalate)

Item No. 35867

CAS Registry No.: 1428862-33-2
Formal Name: 3-[5-[1-(3-methoxypropyl)-4-piperidinyl]-1,3,4-oxadiazol-2-yl]-1-(1-methylethyl)-1H-indazol monoethanedioate
Synonym: SUVN-D4010
MF: $C_{21}H_{29}N_5O_2 \cdot C_2H_2O_4$
FW: 473.5
Purity: $\geq 98\%$
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

Usmarapride (oxalate) is supplied as a solid. A stock solution may be made by dissolving the usmarapride (oxalate) in the solvent of choice, which should be purged with an inert gas. Usmarapride (oxalate) is soluble in methanol and DMSO.

Description

Usmarapride is a partial agonist of the serotonin (5-HT) receptor subtype $5\text{-HT}_{4,1}$. It induces reporter gene expression in CHO cells expressing recombinant human 5-HT_{4E} receptors ($EC_{50} = 44$ nM). *In vivo*, usmarapride (1 and 3 mg/kg) increases novel object exploration time in rats. It also reverses scopolamine-induced memory deficits in rats when administered at doses of 1, 3, and 10 mg/kg.

Reference

1. Nirogi, R., Mohammed, A.R., Shinde, A.K., *et al.* Discovery and preclinical characterization of usmarapride (SUVN-D4010): A potent, selective 5-HT_4 receptor partial agonist for the treatment of cognitive deficits associated with Alzheimer's disease. *J. Med. Chem.* **64**(15), 10641-10665 (2021).

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, 02/22/2024

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