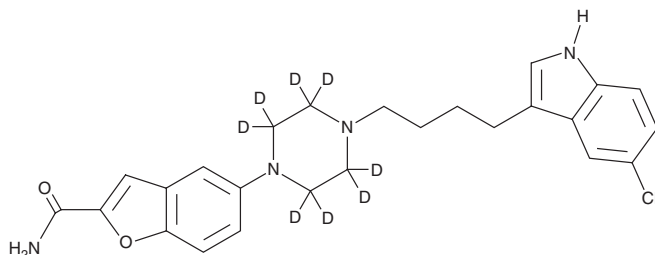


PRODUCT INFORMATION



Vilazodone-d₈ Item No. 25712

CAS Registry No.: 1794789-93-7
Formal Name: 5-(4-(4-(5-cyano-1H-indol-3-yl)butyl)piperazin-1-yl-2,2,3,3,5,5,6,6-d₈)benzofuran-2-carboxamide
MF: C₂₆H₁₉D₈N₅O₂
FW: 449.6
Chemical Purity: ≥98% (Vilazodone)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀
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

Vilazodone-d₈ is intended for use as an internal standard for the quantification of vilazodone (Item No. 21547) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Description

Vilazodone is a selective serotonin reuptake inhibitor (SSRI) and a partial agonist of the serotonin (5-HT) receptor subtype 5-HT_{1A} (IC₅₀s = 0.2 and 0.5 nM, respectively).¹ It increases extracellular 5-HT in the rat ventral hippocampus and frontal cortex when administered intraperitoneally at doses of 1 and 3 mg/kg, respectively. Vilazodone (1 mg/kg, i.p.) decreases immobility in the forced swim test in both rats and mice. Formulations containing vilazodone have been used in the treatment of depression.

Reference

1. Page, M.E., Cryan, J.F., Sullivan, A., *et al.* Behavioral and neurochemical effects of 5-(4-[4-(5-Cyano-3-indolyl)-butyl]-butyl)-1-piperazinyl)-benzofuran-2-carboxamide (EMD 68843): A combined selective inhibitor of serotonin reuptake and 5-hydroxytryptamine_{1A} receptor partial agonist. *J. Pharmacol. Exp. Ther.* **302**(3), 1220-1227 (2002).

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, 12/20/2022

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