

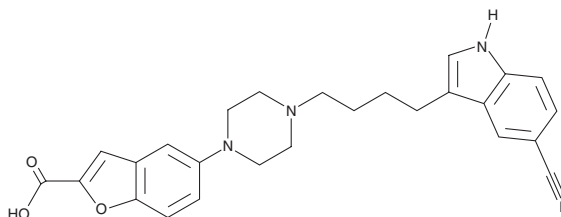
PRODUCT INFORMATION



Vilazodone Carboxylic Acid

Item No. 45834

CAS Registry No.: 163521-19-5
Formal Name: 5-[4-[4-(5-cyano-1H-indol-3-yl)butyl]-1-piperazinyl]-2-benzofurancarboxylic acid
MF: $C_{26}H_{26}N_4O_3$
FW: 442.5
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

Vilazodone carboxylic acid is supplied as a solid. A stock solution may be made by dissolving the vilazodone carboxylic acid in the solvent of choice, which should be purged with an inert gas. Vilazodone carboxylic acid is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Vilazodone carboxylic acid is a metabolite of the selective serotonin reuptake inhibitor (SSRI) vilazodone (Item No. 21547).¹ It is also a potential impurity in commercial preparations of vilazodone.²

References

1. Chavan, B.B., Kalariya, P.D., Tiwari, S., *et al.* Identification and characterization of vilazodone metabolites in rats and microsomes by ultrahigh-performance liquid chromatography/quadrupole time-of-flight tandem mass spectrometry. *Rapid Commun. Mass Spectrom.* **31(23)**, 1974-1984 (2017).
2. Das, P., Srivastava, B., Joseph, S., *et al.* Amidation and N-boc deprotection process improvement for the preparation of 5-(1-piperazinyl)benzofuran-2-carboxamide, a key intermediate of vilazodone. *Org. Process Res. Dev.* **18**, 665-667 (2014).

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, 08/03/2026

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