

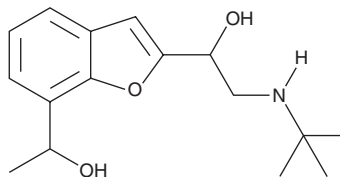
PRODUCT INFORMATION



1'-hydroxy Bufuralol

Item No. 22160

CAS Registry No.: 57704-16-2
Formal Name: α^2 -[[[(1,1-dimethylethyl)amino]methyl]- α^7 -methyl-2,7-benzofurandimethanol
Synonym: Ro 03-7410
MF: $C_{16}H_{23}NO_3$
FW: 277.4
Purity: $\geq 97\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Special Conditions: Hygroscopic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

1'-hydroxy Bufuralol is supplied as a solid. A stock solution may be made by dissolving the 1'-hydroxy bufuralol in the solvent of choice. 1'-hydroxy Bufuralol is soluble in the organic solvent chloroform, which should be purged with an inert gas. 1'-hydroxy Bufuralol is slightly soluble in methanol.

Description

1'-hydroxy Bufuralol is a metabolite of bufuralol (Item No. 17794). Bufuralol is metabolized by the cytochrome P450 isoforms 2D1 and 2D2 (CYP2D1/2); therefore, quantification of the bufuralol metabolite, 1'-hydroxy bufuralol, can be used to determine CYP2D1/2 enzyme activity in humans and rats.^{1,2}

References

1. Daskalopoulos, E.P., Lang, M.A., Marselos, M., *et al.* D₂-dopaminergic receptor-linked pathways: Critical regulators of CYP3A, CYP2C, and CYP2D. *Mol. Pharmacol.* **82**(4), 668-678 (2012).
2. Yamazaki, H., Guo, Z., Persmark, M., *et al.* Bufuralol hydroxylation by cytochrome P450 2D6 and 1A2 enzymes in human liver microsomes. *Mol. Pharmacol.* **46**(3), 568-577 (1994).

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, 12/07/2022

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