

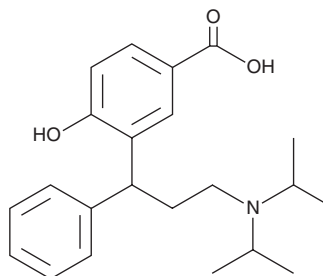
# PRODUCT INFORMATION



## *rac*-5-carboxy Tolterodine

Item No. 34074

**CAS Registry No.:** 1076199-77-3  
**Formal Name:** 3-[3-[*bis*(1-methylethyl)amino]-1-phenylpropyl]-4-hydroxy-benzoic acid  
**MF:** C<sub>22</sub>H<sub>29</sub>NO<sub>3</sub>  
**FW:** 355.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 253 nm  
**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

*rac*-5-carboxy Tolterodine is supplied as a solid. A stock solution may be made by dissolving the *rac*-5-carboxy tolterodine in the solvent of choice, which should be purged with an inert gas. *rac*-5-carboxy Tolterodine is soluble in ethanol.

### Description

*rac*-5-carboxy Tolterodine is an inactive metabolite of the muscarinic acetylcholine receptor antagonist tolterodine (Item No. 15027).<sup>1</sup> It is formed from tolterodine via a 5-hydroxymethyl tolterodine (Item No. 27833) intermediate.

### Reference

1. Clemett, D. and Jarvis, B. Tolterodine: A review of its use in the treatment of overactive bladder. *Drugs Aging* **18**(4), 277-304 (2001).

#### 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, 11/17/2022

#### 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