

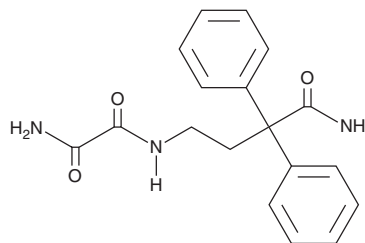
PRODUCT INFORMATION



Imidafenacin Metabolite M4

Item No. 34301

CAS Registry No.: 503598-17-2
Formal Name: N¹-(4-amino-4-oxo-3,3-diphenylbutyl)-ethanediamide
MF: C₁₈H₁₉N₃O₃
FW: 325.4
Purity: ≥90%
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

Imidafenacin metabolite M4 is supplied as a solid. A stock solution may be made by dissolving the imidafenacin metabolite M4 in the solvent of choice, which should be purged with an inert gas. Imidafenacin metabolite M4 is soluble in the organic solvent methanol.

Description

Imidafenacin metabolite M4 is a metabolite of the muscarinic acetylcholine receptor antagonist imidafenacin.¹ It is formed from imidafenacin by the cytochrome P450 (CYP) isoform CYP3A4.

Reference

1. Kanayama, N., Kanari, C., Masuda, Y., *et al.* Drug-drug interactions in the metabolism of imidafenacin: Role of the human cytochrome P450 enzymes and UDP-glucuronic acid transferases, and potential of imidafenacin to inhibit human cytochrome P450 enzymes. *Xenobiotica* **37(2)**, 139-154 (2007).

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, 11/18/2022

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