

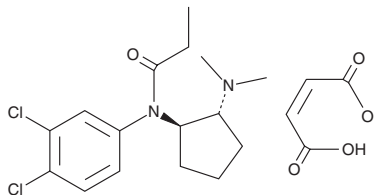
PRODUCT INFORMATION



U-48753E (maleate)

Item No. 20734

CAS Registry No.: 67450-45-7
Formal Name: *trans*-N-(3,4-dichlorophenyl)-N-[(1*R*,2*R*)-2-(dimethylamino)cyclopentyl]-propanamide, 2*Z*-butenedioate
MF: C₁₆H₂₂Cl₂N₂O • C₄H₄O₄
FW: 445.3
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

U-48753E (maleate) (Item No. 20734) is an analytical reference standard that is functionally categorized as a nontricyclic antidepressant. It has been shown to modulate signaling through yohimbine (Item No. 19869), oxotremorine, and apomorphine in mice, although its physiological and toxicological effects in humans are not known.¹ This product is intended for research and forensic applications.

Reference

1. Szmuszkovicz, J., Von Voigtlander, P.F., and Kane, M.P. A new nontricyclic antidepressant agent. Synthesis and activity of N-[*trans*-2-(dimethylamino)cyclopentyl]-N-(3,4-dichlorophenyl)propanamide and related compounds. *J. Med. Chem.* **24**(10), 1230-1236 (1981).

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, 06/22/2023

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