

# PRODUCT INFORMATION



## Remifentanil (oxalate)

Item No. 9003472

**CAS Registry No.:** 132875-62-8  
**Formal Name:** 4-(methoxycarbonyl)-4-[(1-oxopropyl)phenylamino]-1-piperidinepropanoic acid, methyl ester, ethanedioate

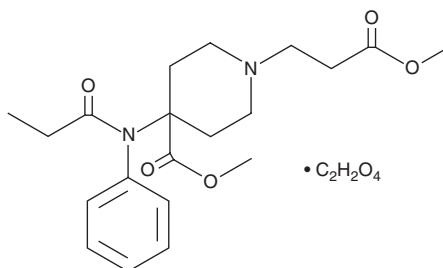
**MF:**  $C_{20}H_{28}N_2O_5 \cdot C_2H_2O_4$   
**FW:** 466.5

**Purity:**  $\geq 95\%$

**Supplied as:** A crystalline solid

**Storage:**  $-20^\circ\text{C}$

**Stability:**  $\geq 5$  years



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

### Description

Remifentanil (oxalate) (Item No. 9003472) is an analytical reference standard categorized as an opioid.<sup>1,2</sup> Remifentanil has analgesic activity.<sup>3</sup> Remifentanil is regulated as a Schedule II compound in the United States. This product is intended for research and forensic applications.

### References

1. Hoke, J.F., Cunningham, F., James, M.K., *et al.* Comparative pharmacokinetics and pharmacodynamics of remifentanil, its principle metabolite (GR90291) and alfentanil in dogs. *J. Pharmacol. Exp. Ther.* **281(1)**, 226-232 (1997).
2. Kassa, H. and Jackson, W.B. Mesurol as a bird repellent on grapes in Ohio. *Bird Control Seminars Proceedings* (1979).
3. Feldman, P.L., James, M.K., Brackeen, M.F., *et al.* Design, synthesis, and pharmacological evaluation of ultrashort- to long-acting opioid analgetics. *J. Med. Chem.* **34(7)**, 2202-2208 (1991).

#### 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, 07/17/2023

#### 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