

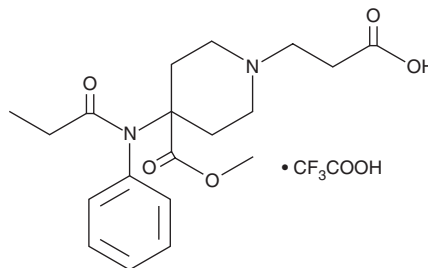
PRODUCT INFORMATION



Remifentanil Acid (trifluoroacetate salt)

Item No. 21926

CAS Registry No.: 132875-69-5
Formal Name: 4-(methoxycarbonyl)-4-[(1-oxopropyl)phenylamino]-1-piperidinepropanoic acid, 2,2,2-trifluoroacetate
Synonyms: GI-90291, GR 90291
MF: C₁₉H₂₆N₂O₅ • CF₃COOH
FW: 476.4
Purity: ≥98%
UV/Vis.: λ_{max}: 205 nm
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

Remifentanil acid (trifluoroacetate salt) (Item No. 21926) is an analytical reference standard categorized as an opioid.¹ It is a known active metabolite of remifentanil (Item No. 19291).² This product is intended for research and forensic applications.

Reference

1. 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).
2. 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).

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, 07/06/2023

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