

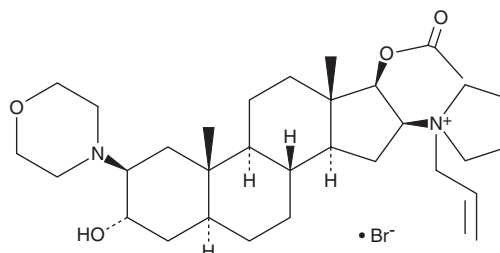
PRODUCT INFORMATION



Rocuronium (bromide)

Item No. 23698

CAS Registry No.: 119302-91-9
Formal Name: 1-[(2 β ,3 α ,5 α ,16 β ,17 β)-17-(acetyloxy)-3-hydroxy-2-(4-morpholinyl)androstan-16-yl]-1-(2-propenyl)-pyrrolidinium, monobromide
Synonym: (+)-Rocuronium
MF: C₃₂H₅₃N₂O₄ • Br
FW: 609.7
Purity: ≥95%
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

Rocuronium (bromide) is supplied as a solid. A stock solution may be made by dissolving the rocuronium (bromide) in the solvent of choice. Rocuronium (bromide) is soluble in the organic solvent chloroform, which should be purged with an inert gas. It is also soluble in water.

Description

Rocuronium is a non-depolarizing muscle relaxant.¹ It selectively induces neuromuscular blockade in cats (ED₅₀s = 153, 185, 1,329, and 4,119 µg/kg for tibialis, soleus, and vagus muscles, and nictitating membranes, respectively) with a more rapid onset time than vecuronium (Item No. 15603). Formulations containing rocuronium have been used for surgical anesthesia.²

References

1. Muir, A.W., Houston, J., Green, K.L., *et al.* Effects of a new neuromuscular blocking agent (Org 9426) in anaesthetized cats and pigs and in isolated nerve-muscle preparations. *Br. J. Anaesth.* **63**(4), 400-410 (1989).
2. Fawcett, W.J., Fauvel, N.J., and Feldman, S.A. Comparison of recovery index of rocuronium or vecuronium with simultaneously administered pancuronium in the isolated forearm. *Anaesthesia* **48**(3), 200-201 (1993).

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, 04/03/2024

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