

PRODUCT INFORMATION

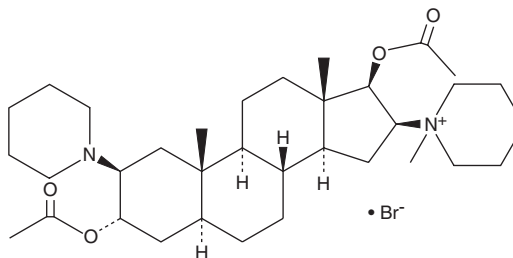


Vecuronium (bromide)

Item No. 15603

CAS Registry No.: 50700-72-6
Formal Name: 1-[(2 β ,3 α ,5 α ,16 β ,17 β)-3,17-bis(acetyloxy)-2-(1-piperidinyl)androstan-16-yl]-1-methylpiperidinium, monobromide

Synonym: NC 45
MF: C₃₄H₅₇N₂O₄ • Br
FW: 637.7
Purity: ≥98%
Supplied as: A solid
Storage: Room temperature
Stability: ≥4 years



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

Laboratory Procedures

Vecuronium (bromide) is supplied as a solid. A stock solution may be made by dissolving the vecuronium (bromide) in the solvent of choice. Vecuronium (bromide) is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas. The solubility of vecuronium (bromide) in these solvents is approximately 100 mM. Vecuronium (bromide) is also soluble in water at a concentration of approximately 25 mM.

Description

Vecuronium is a non-depolarizing muscle relaxant derived from the aminosteroid pancuronium (Item No. 23778) and used adjunctively to general anesthesia.¹ It competitively blocks cholinergic receptors at the motor end plate of the neuromuscular junction, inducing temporary paralysis.² In humans, it has been shown to reduce muscle twitch tension with an ED₅₀ value of 0.15 mg/kg for a duration of 27 minutes without inducing cardiovascular effects.¹

References

1. Fahey, M.R., Morris, R.B., Miller, R.D., *et al.* Clinical pharmacology of ORG NC45 (Norcuron™): A new nondepolarizing muscle relaxant. *Anesthesiology* **55**(1), 6-11 (1981).
2. Raghavendra, T. Neuromuscular blocking drugs: Discovery and development. *J. R. Soc. Med.* **95**(7), 363-367 (2002).

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, 10/03/2018

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