

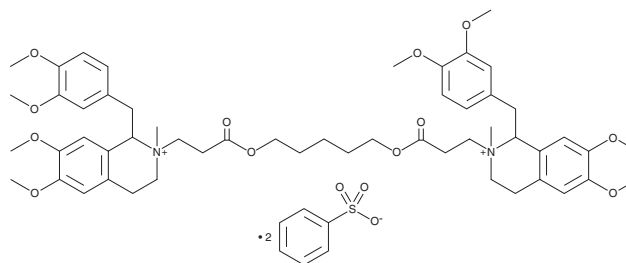
PRODUCT INFORMATION



Atracurium (besylate)

Item No. 17796

CAS Registry No.: 64228-81-5
Formal Name: 2,2'-[1,5-pentanediy]bis[oxy(3-oxo-3,1-propanediy)]bis[1-[(3,4-dimethoxyphenyl)methyl]-1,2,3,4-tetrahydro-6,7-dimethoxy-2-methyl-isoquinolinium, dibenzenesulfonate
Synonyms: BW 33A, Tracrium, Wellcome 33A74
MF: C₅₃H₇₂N₂O₁₂ • 2C₆H₅O₃S
FW: 1,243.5
Purity: ≥95%
UV/Vis.: λ_{max}: 282 nm
Supplied as: A crystalline 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

Atracurium (besylate) is supplied as a crystalline solid. A stock solution may be made by dissolving the atracurium (besylate) in the solvent of choice. Atracurium (besylate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of atracurium (besylate) in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of atracurium (besylate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of atracurium (besylate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Atracurium (besylate) is a benzylisoquinoline that acts as a neuromuscular blocking agent with slow onset and intermediate duration.^{1,2} It causes non-depolarizing block of the neuromuscular junction.²

References

1. Mirakhur, R. New developments in nondepolarizing muscle relaxants. *Yale J. Biol. Med.* **66**, 463-471 (1994).
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, 01/05/2023

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