

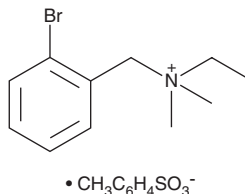
PRODUCT INFORMATION



Bretylum (tosylate)

Item No. 29979

CAS Registry No.: 61-75-6
Formal Name: 2-bromo-N-ethyl-N,N-dimethylbenzenemethanaminium, 4-methylbenzenesulfonate (1:1)
MF: C₁₁H₁₇BrN • C₇H₇O₃S
FW: 414.4
Purity: ≥95%
UV/Vis.: λ_{max}: 220 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

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

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 bretylum (tosylate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of bretylum (tosylate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Bretylum is a class III antiarrhythmic agent and an inhibitor of the Na⁺/K⁺-ATPase (IC₅₀ = 4.5 mM).¹ Bretylum also has antiadrenergic activity, inhibiting auricular nerve stimulation-induced vasoconstriction in isolated rabbit ears and hypogastric nerve stimulation-induced contraction of isolated rabbit uterus.² It inhibits neuroeffector calcium transients (NCTs), as well as increases action potential delay and the absolute refractory period, but does not inhibit field stimulus-induced CTs in isolated mouse vas deferens sympathetic nerve terminals.³ Bretylum prevents ventricular fibrillation in anesthetized dogs in a model of sudden coronary death when administered at a dose of 10 mg/kg.⁴ Formulations containing bretylum were previously used in the prevention and treatment of ventricular fibrillation.

References

1. Helms, J.B., Arnett, K.L., Gatto, C., *et al.* Bretylum, an organic quaternary amine, inhibits the Na,K-ATPase by binding to the extracellular K-site. *Blood Cells Mol. Dis.* **32(3)**, 394-400 (2004).
2. Boura, A.L., Copp, F.C., and Green, A.F. New antiadrenergic compounds. *Nature* **184**, BA70-BA71 (1959).
3. Brain, K.L. and Cunnane, T.C. Bretylum abolishes neurotransmitter release without necessarily abolishing the nerve terminal action potential in sympathetic terminals. *Br. J. Pharmacol.* **153(4)**, 831-839 (2007).
4. Holland, K., Patterson, E., and Lucchesi, B.R. Prevention of ventricular fibrillation by bretylum in a conscious canine model of sudden coronary death. *Am. Heart J.* **105(5)**, 711-717 (1983).

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, 05/06/2020

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