

PRODUCT DATA SHEET



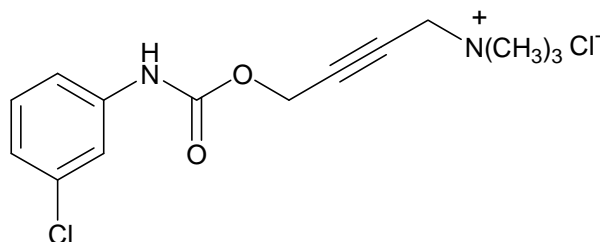
PRODUCT: McN-A-343
4-(N-[3-Chlorophenyl]carbamoxyloxy)-2-butynyl-
trimethylammonium chloride

CAS NO.: 55-45-8

CATALOG NO.: C-119

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: $C_{14}H_{18}ClN_2O_2 \cdot Cl$

MOLECULAR WEIGHT: 317.2

PURITY: >98%

SOLUBILITY: Soluble in water (>25 mg/mL)

STORAGE: Store, as supplied, at room temperature for up to 1 year. Store solutions at $-20^{\circ}C$ for up to 3 months.

APPLICATION NOTES: A potent and selective muscarinic M1 receptor agonist^{1,2}. $10^{-3} - 10^{-5}$ M concentrations have been shown to induce receptor response in mouse stomach fundus³. This compound has also been shown to have opposing effects on the M3 receptor, which in some situations may mask M1 receptor response³.

REFERENCES:

- 1) R.H. Davies *et al.* *J. Pharm. Pharmacol.* 2001 **53** 487
- 2) J. Liu *et al.* *J. Pharmacol. Exp. Ther.* 2002 **302** 397
- 3) P.W. Stengel and M.L. Cohen *J. Pharmacol. Exp. Ther.* 2003 **304** 675

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.