

PRODUCT DATA SHEET



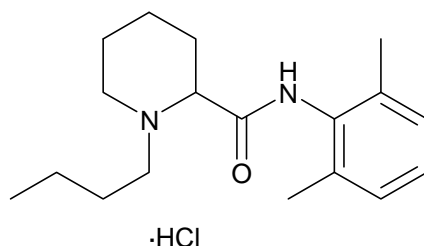
PRODUCT: Bupivacaine HCl
1-Butyl-N-(2,6-dimethylphenyl)-2-piperidine
carboxamide·HCl

CAS NO.: 14252-80-3

CATALOG NO.: NA-139

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: $C_{18}H_{28}N_2O \cdot HCl$

MOLECULAR WEIGHT: 324.9

PURITY: 99% (TLC: 1% Methanol/methylene chloride + 0.5% NH_4OH ; $R_f = 0.35$)

SOLUBILITY: Soluble in: Water (5 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: Sodium channel blocker¹. In rat dorsal horn neurons externally applied bupivacaine produced tonic block of Na^+ current with an $IC_{50}=26 \mu M^1$. Clinically useful local anesthetic². Also inhibits TASK ($IC_{50}=68 \mu M^3$)

REFERENCES: 1. A. Olschewski *et al. Anesthesiology* 1998 **88** 172
2. H. Bariskaner *et al. Methods Find. Exp. Clin. Pharmacol.* 2007 **29** 337
3. D. Leonoudakis *et al. J. Neurosci.* 1998 **18** 868

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.