

PRODUCT DATA SHEET



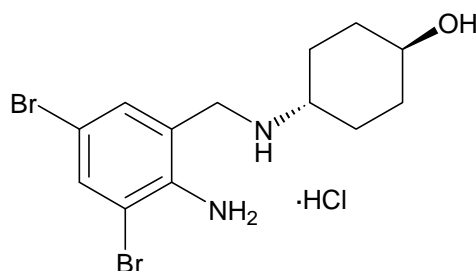
PRODUCT: Ambroxol
trans-4-(2-Amino-3,5-dibromobenzylamino)
cyclohexane·HCl

CAS NO. 18683-91-5

CATALOG NO.: FR-112

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: C₁₃H₁₈Br₂N₂O·HCl

MOLECULAR WEIGHT: 414.6

PURITY: 98%(TLC)

SOLUBILITY: Soluble in water

STORAGE: Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.

APPLICATION NOTES: Radical scavenger. Inhibits doxorubicin-induced lipid peroxidation in vivo (iv 70 mg/kg, rat).¹ Blocks generation of reactive oxygen species by bronchoalveolar lavage cells.² Anti-inflammatory activity: inhibits the release of inflammatory mediators (@100 μM) from human leukocytes and mast cells.³

REFERENCES: 1. D.Nowak *et al. Free Radic.Biol.Med.* 1995 **19** 659
2. S.Teramoto *et al. Pharmacology* 1999 **59** 135
3. B.F.Gibbs *et al. Inflamm.Res.* 1999 **48** 86

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.