

PRODUCT DATA SHEET



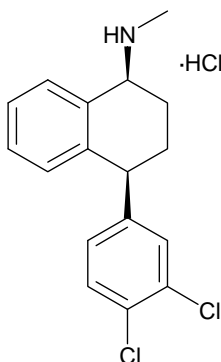
PRODUCT: Sertraline HCl
CP-51974
(1S,4S)-4-(3,4-Dichlorophenyl)-1,2,3,4-tetrahydro-N-methyl-1-naphthalenamine hydrochloride

CAS NO.: 79559-97-0

CATALOG NO.: NS-115

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: C₁₇H₁₇Cl₂N·HCl

MOLECULAR WEIGHT: 342.7

PURITY: 98% (TLC)

SOLUBILITY: Soluble in: DMSO (>25 mg/mL); ethanol (10 mg/mL); water (3 mg/mL)

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

APPLICATION NOTES: A selective serotonin transporter inhibitor¹. In clinical use as an antidepressant². Displays antiproliferative activity on human colorectal carcinoma cells³.

REFERENCES: 1. B.K. Koe. *J. Pharmacol. Exp. Ther.* 1983 **226** 686
2. H.R. Khouzam *et al. Compr. Ther.* 2003 **29** 47
3. I. Gil-ad *et al. Int. J. Oncol.* 2008 **33** 277

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.