

PRODUCT DATA SHEET



PRODUCT: ORG-12962
1-[6-Chloro-5-trifluoromethyl]-2-pyridinyl]-
piperazine

CAS NO.: 210821-63-9

CATALOG NO.: NS-705

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: Pale yellow solid

MOLECULAR FORMULA: C₁₀H₁₁ClF₃N₃

MOLECULAR WEIGHT: 265.7

PURITY: 98% (TLC: 5% MeOH/methylene chloride+1% NH₄OH; R_f = 0.33)

SOLUBILITY: Soluble in: DMSO (>25 mg/mL); Ethanol (>25 mg/mL)

APPLICATION NOTES: ORG-12962 is a selective serotonin 5-HT_{2C} agonist¹. It displays antiaversive effects in a panic-like anxiety rat model². May display appetite regulating properties similar to other 5-HT_{2C} agonists³.

REFERENCES: 1. R.H. Porter *et al. Br. J. Pharmacol.* 1999 **128** 13
2. F. Jenck *et al. Eur. Neuropsychopharmacol.* 1998 **8** 161
3. J.C. Halford *et al. Curr. Drug Targets* 2005 **6** 201

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

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.