

PRODUCT DATA SHEET



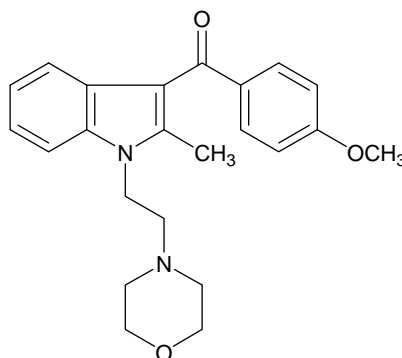
PRODUCT: Pravadoline
(4-Methoxyphenyl)(2-methyl-1-(2-(4-morpholinyl)ethyl)-1H-indol-3-yl)methanone

CAS NO: 92623-83-1

CATALOG NO.: CR-110

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: Off-white solid (m.p. 107-110°)

MOLECULAR FORMULA: $C_{23}H_{26}N_2O_3$

MOLECULAR WEIGHT: 378.5

PURITY: 98%

SOLUBILITY: Soluble in DMSO

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

APPLICATION NOTES: Pravadoline is a novel analgesic agent whose mechanism of action includes inhibition of cyclooxygenase¹ and cannabinoid receptor agonist activity². The product is not sterile.

REFERENCES: 1. D.R.Haubrich *et al. J.Pharmacol.Exp.Ther.* 1990 **255** 511
2. T.E.D'Ambra *et al. J.Med.Chem.* 1992 **35** 124

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.