

PRODUCT DATA SHEET



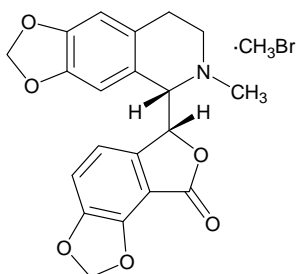
PRODUCT: (-)-Bicuculline methobromide

CAS NO: 73604-30-5

CATALOG NO.: EA-109

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: $C_{21}H_{20}BrNO_6$

MOLECULAR WEIGHT: 462.3

PURITY: 98%

SOLUBILITY: Soluble in water (25 mg/ml)

STORAGE: Store, as supplied, at 0-4°C for up to 1 year. Store solutions at -20°C for up to 3 months.

APPLICATION NOTES: GABA_A antagonist.¹

REFERENCES: 1. D.R.Hill and N.G.Bowery *Nature* 1981 **290** 149

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.