

PRODUCT DATA SHEET



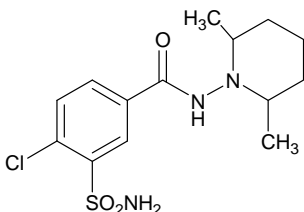
PRODUCT: Clopamide
4-Chloro-N-(2,6-dimethylpiperidinyl)-
3-sulfamoylbenzamide

CAS NO: 636-54-4

CATALOG NO.: CM-112

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 248-249°C)

MOLECULAR FORMULA: $C_{14}H_{20}ClN_3O_3S$ MOLECULAR WEIGHT: 345.8

PURITY: 98%(TLC)

SOLUBILITY: Soluble in DMSO (20 mg/mL) and ethanol (5 mg/mL warm)

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

APPLICATION NOTES: Clinically useful diuretic.¹ The product is not sterile.

REFERENCES: 1. J.J.McNeil *et al. Clin.Pharmacol.Ther.* 1987 **42** 299

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.