

PRODUCT DATA SHEET



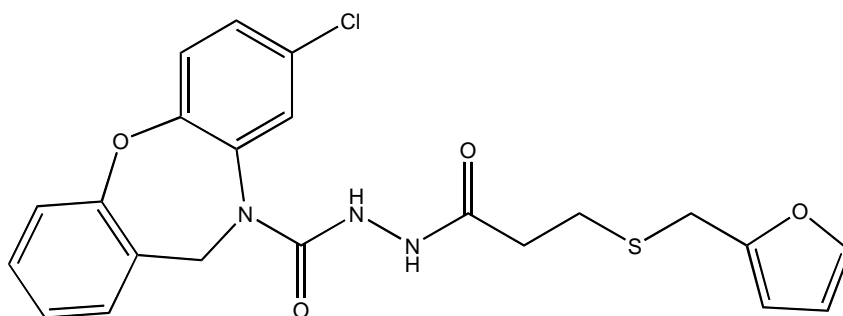
PRODUCT: SC-51322

CAS.NO:

CATALOG NO.: RA-112

LOT#: temp

STRUCTURE:



PHYSICAL APPEARANCE: white solid

MOLECULAR FORMULA: $C_{22}H_{20}ClN_3O_4S$

MOLECULAR WEIGHT: 457.9

PURITY: 98% (TLC)

SOLUBILITY: Soluble in ethanol (10mg/ml), DMSO (25mg/ml)

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

APPLICATION NOTES: Prostaglandin E₂ (EP1 receptor) antagonist (pA₂=8.1 in guinea pig ileum muscle strip assay)^{1,2}. Has *in vivo* analgesic activity (ED₅₀=0.9 mg/kg, mouse)^{1,2}. Does not inhibit cyclooxygenase².

REFERENCES: 1. Hallinan, EA *et al. J. Med. Chem.* 1996 **39** 609
2. E.A. Hallinan *et al. Bioorg. Med. Chem. Lett.* 1994 **4** 509

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.