

PRODUCT DATA SHEET



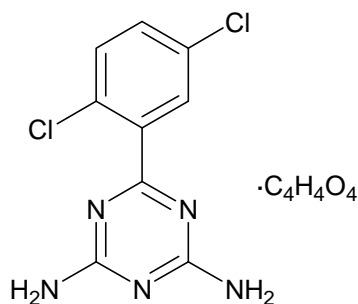
PRODUCT: Irsogladine maleate

CAS NO. 84504-69-8

CATALOG NO.: PD-141

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: $C_9H_7Cl_2N_5 \cdot C_4H_4O_4$ MOLECULAR WEIGHT: 372.2

PURITY: 98% (TLC: 10% Methanol/methylene chloride + 0.5% NH₄OH; R_f = 0.45)

SOLUBILITY: Soluble in DMSO (25 mg/mL)

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

APPLICATION NOTES: Irsogladine is a selective PDE4 inhibitor¹ which also enhances gap junction intercellular communication² via upregulation of connexins 32 and 263. It inhibits angiogenesis in wild-type and plasminogen activator-deficient mice.⁴

- REFERENCES:
1. T. Kyojima *et al.* *Life Sci.* 2004 **75** 1833
 2. Y. Kawasaki *et al.* *Pancreas* 2002 **25** 373
 3. Y. Nakashima *et al.* *Hepatol. Res.* 2000 **18** 29
 4. C. J. Ren *et al.* *J. Surg. Res.* 1998 **77** 126

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.