

PRODUCT DATA SHEET



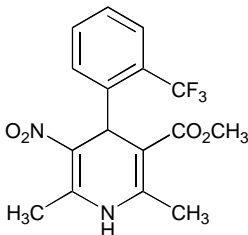
PRODUCT: BAY-K8644

CAS NO: 98791-67-4

CATALOG NO.: CA-410

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: Yellow solid (m.p. 168-170°C)

MOLECULAR FORMULA: C₁₆H₁₅F₃N₂O₄

MOLECULAR WEIGHT: 356.3

PURITY: 98%

SOLUBILITY: Soluble in ethanol (20 mg/ml) and DMSO (20 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: Potent, direct acting voltage-sensitive Ca²⁺ channel activator.^{1,2} It exerts positive inotropic, vasoconstrictive and positive chronotropic activity (*in vitro* and *in vivo*) in a concentration range from 10⁻⁹ to 10⁻⁷ M³.

REFERENCES: 1. J.G.Sarmiento *et al.* *J.Pharmacol.Exp.Ther.* 1987 **241** 140
2. G.Thomas *et al.* *Circulation Res.* 1985 **56** 87
3. P.L.Vaghy *et al.* *Circulation Res.* 1984 **55** 549

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.