

# PRODUCT DATA SHEET



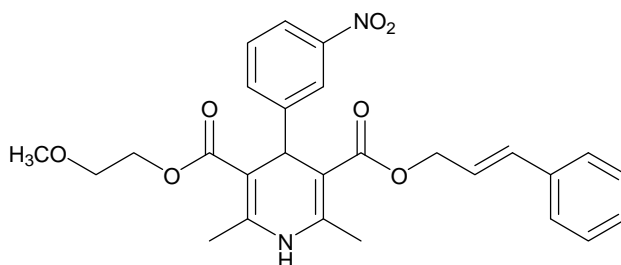
PRODUCT: Cilnidipine  
FRC-8653  
1,4-Dihydro-2,6-dimethyl-4-(3-nitrophenyl)-3,5-pyridinedicarboxylic acid 2-methoxyethyl (2E)-3-phenyl-2-propenyl ester

CAS NO.: 132203-70-4

CATALOG NO.: CA-237

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA:  $C_{27}H_{28}N_2O_7$

MOLECULAR WEIGHT: 492.5

PURITY: 98% (TLC: 5% Methanol/methylene chloride;  $R_f = 0.54$ )

SOLUBILITY: Soluble in: DMSO (>25 mg/mL); ethanol (20 mg/mL)

STORAGE: Store as supplied, at room temperature for up to 1 year. Store solutions at  $-20^{\circ}\text{C}$  for up to 3 months.

APPLICATION NOTES: Dual L- / N-type calcium channel blocker<sup>1</sup>. Displays renal and vascular protective effects in patients with essential hypertension<sup>2</sup>. Displays neuroprotective effects in a rat focal brain ischemia model<sup>3</sup>.

REFERENCES: 1. K. Hayashi et al. *Circ. Res.* 2007 **100** 342  
2. S. Morimoto et al. *J. Hypertens.* 2007 **25** 2178  
3. A. Takahara et al. *Biol. Pharm. Bull.* 2004 **27** 1388

**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.**