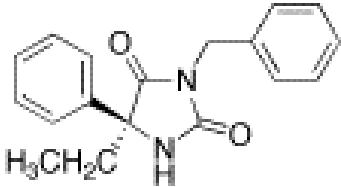


Potent and selective inhibitor of CYP2C19

## Instruction Manual

Catalog Number	PK-CA577-2920-5
Description	It is a potent and selective CYP2C19 inhibitor. Our product is the S enantiomer and not a racemic mixture.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	294.35 g/mol; C <sub>18</sub> H <sub>18</sub> N <sub>2</sub> O <sub>2</sub>
Chemical Structure	
Purity	≥98%
Appearance / Formulation / Solubility	Off-white solid; DMSO (>20 mg/ml)
Storage & Stability	Stable at -20 °C for 2 years
Applications	See description
References	Drug Metab. Dispos., 30, 235 (2002), Drug Metab. Dispos., 31, 343 (2003)
Caution	Protect from air, light and moisture. Do not take internally. Wear gloves and mask when handling since this product is toxic to the skin or eyes!

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.