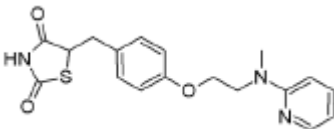


Instruction Manual

Catalog Number	PK-CA577-1559-100
Description	Belongs to the thiazolidinedione class of anti-diabetic drugs. A selective ligand of PPAR γ that has no PPAR α -binding action. Induces apoptosis in various cells, e.g. in vascular smooth muscle cells.
Quantity	100 mg
Sequence / Molecular Weight / Molecular Formula	357.4 Da; C ₁₅ H ₁₉ N ₃ O ₃ S
Purity	>99%
Appearance / Formulation / Solubility	White to off-white powder. Soluble in DMSO (>30 mg/ml).
Chemical Structure	 The chemical structure of Rosiglitazone consists of a thiazolidinedione ring system (a five-membered ring with a sulfur atom and two carbonyl groups, one of which is attached to an NH group). This ring is connected via a methylene group to a para-substituted benzene ring. The benzene ring is further connected via an ether linkage to a propyl chain, which is terminated by a dimethylamino group (-N(CH ₃) ₂). The dimethylamino group is attached to a pyridine ring.
Storage & Stability	Store at -20°C in a dry environment.
Applications	see Description
References	NA
Caution	Do not ingest, swallow, or inhale. Do not get in eyes or on skin. Care should be taken to prevent contact through all routes of exposure. Toxicity information is not available at this time.

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