

Tyrphostin SU 1498; (E)-N-(3''-Phenylpropyl)- α -cyano-3',5'-diisopropyl-4'-hydroxycinnamamide

Instruction Manual

Catalog Number	PK-CA577-1836-5
Description	A potent and selective inhibitor (IC ₅₀ = 700 nM) of vascular endothelial growth factor receptor 2 (VEGFR 2) (also known as Flk1). Very weak inhibitor of PDGFR-kinase (IC ₅₀ >50 μ M), EGFR-kinase and HER-2 kinase. SU1498 has been shown to cause accumulation of phosphorylated ERK and inhibits its activity in vivo and in vitro by inhibition of ERK dephosphorylation.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	390.52 Da; C ₂₅ H ₃₀ N ₂ O ₂
Chemical Structure	
Purity	≥ 98% as determined by TLC
Appearance / Formulation / Solubility	Yellow Solid; Soluble in DMSO (~20 mg/ml)
Storage & Stability	Store at -20°C.
Applications	see Description
References	Boguslawski, G., et al. (2004). J. Biol. Chem. 279, 5716-5724.
Caution	Do not take internally. Wear gloves and mask when handling the product! Avoid contact by all modes of exposure.

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