

(3Z)-N-(3-Chlorophenyl)-3-({3,5-dimethyl-4-[(4-methylpiperazin-1-yl)carbonyl]-1H-pyrrol-2-yl}methylene)-N-methyl-2-oxo-2,3-dihydro-1H-indole-5-sulfonamide**Instruction Manual**

Catalog Number	PK-CA577-1938-25
Description	A specific, ATP-competitive small molecule inhibitor of the catalytic activity of Met. Displays selectivity for Met enzyme versus a panel of other tyrosine kinases with the following IC50 values: Met = 0.02 μ M, Flk = 1.3 μ M, EGFR = >100 μ M, PDGF β R = >20 μ M, Tie2 = >100 μ M, c-src = >10 μ M, cdk2 = >10 μ M, and FGFR-1 = 9.7 μ M). Inhibition of the Met kinase activity by SU11274 leads to time- and dose-dependent reduced cell growth and induced G1 cell cycle arrest and apoptosis.
Quantity	25 mg
Sequence / Molecular Weight / Molecular Formula	568.09 Da; C ₂₈ H ₃₀ ClN ₅ O ₄ S
Chemical Structure	
Purity	≥98% as determined by HPLC
Appearance / Formulation / Solubility	Orange Solid; Soluble in DMSO (10 mg/ml on warming)
Storage & Stability	Store at -20°C. Protect from light.
Applications	see Description
References	Berthou, S., et al. (2004). <i>Oncogene</i> 23, 5387-5393. Sattler, M., et al. (2003). <i>Cancer Res.</i> 63, 5462-5469.
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.