

Instruction Manual

Catalog Number	PK-CA577-1570-5
Description	Specific and potent inhibitor of JAK2 protein tyrosine kinase, EGF receptor autophosphorylation, DNA synthesis and cell growth. Selectively inhibits EGF receptor autophosphorylation (IC50 = 100 µM) and selectively inhibits DNA synthesis and cell growth; induces apoptosis. Blocks growth of leukemia cells in vivo and in vitro.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	294,31 Da; C ₁₅ H ₁₄ N ₂ O ₃
Purity	>99%
Appearance / Formulation / Solubility	Off-white solid. Soluble in ethanol (5 mg/ml) and DMSO. Insoluble in water.
Storage & Stability	Store at -20°C. Do not expose to sunlight or fluorescent lights for extended periods.
Applications	see Description
References	Abe, et al., J. Biol. Chem. 274, 21003-21010, (1999). Al-Shami, et al., J. Biol. Chem. 274, 5333, (1999). Eriksen, K.W., et al., Leukemia 15, 787-793, (2001). Potier, et al., Brain Res. 816, 587-597, (1999).
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.

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