

2-(4-(4-Chlorophenoxy)phenyl)-1H-benzimidazole-5-carboxamide

Instruction Manual

Catalog Number	PK-CA577-1702
Description	A cell-permeable, potent, and ATP-competitive inhibitor of Chk2 (checkpoint kinase 2; $IC_{50} = 15$ nM; $K_i = 37$ nM). Displays ~1,000-fold greater selectivity for Chk2 compared to Cdk1/cyclin B and CK1 ($IC_{50} = 12$ μ M and 17 μ M, respectively) and only weakly affects the activities of a panel of 31 kinases (< 25% inhibition at 10 μ M), including Chk1. Shown to rescue both peripheral CD4+ and CD8+ T-cells from γ -irradiation induced apoptosis with an EC_{50} of 3 μ M and 7.6 μ M, respectively.
Quantity	1 mg
Sequence / Molecular Weight / Molecular Formula	363.8 Da; $C_{20}H_{14}ClN_3O_2$
Purity	>98% by HPLC analysis
Appearance / Formulation / Solubility	Pale yellow solid. Soluble in DMSO (10 mg/ml).
Storage & Stability	Store at -20°C. Protect from light and moisture. Stable for 1 year as supplied under proper storage conditions.
Applications	see Description
References	NA
Caution	NA

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