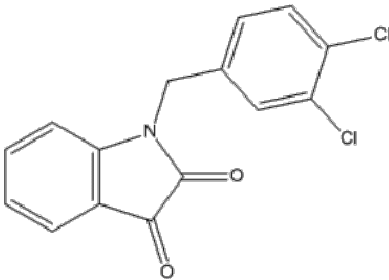


Instruction Manual

Catalog Number	PK-CA577-2784-5
Description	Apoptosis activator 2 is a cell-permeable apoptosis activator that promotes the cytochrome c-dependent oligomerization of Apaf-1 into the mature apoptosome. Increases procaspase-9 processing and subsequent caspase-3 activation. Induces apoptosis in tumor cells ($IC_{50} = 4 - 9 \mu M$ for leukemia cells) with weak or no effect on normal cell lines or those defective/deficient in Apaf-1, caspase-9 or caspase-3 activity ($IC_{50} > 40 \mu M$).
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	306.14 Da; $C_{15}H_9Cl_2NO_2$
Chemical Structure	
Synonyms	1-[(3,4-dichlorophenyl)methyl]indole-2,3-dione
Purity	$\geq 98\%$ as determined by HPLC.
Appearance / Formulation / Solubility	Orange solid. Soluble in DMSO (~30 mg/ml).
Storage & Stability	Store desiccated at $-20^{\circ}C$. Protect from air and light.
Applications	Apoptosis Inducer. See Description.
References	Nguyen, J.T., and Wells, J.A. (2003). Proc. Natl. Acad. Sci. USA 100, 7533-7538.
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.