

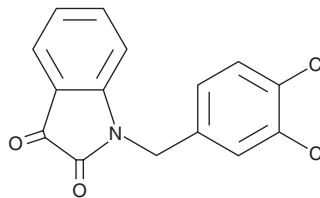
# PRODUCT INFORMATION



## Apoptosis Activator 2

Item No. 10004176

**CAS Registry No.:** 79183-19-0  
**Formal Name:** 1-[(3,4-dichlorophenyl)methyl]-1H-indole-2,3-dione  
**Synonyms:** AAIL, N-(3,4-dichlorobenzyl) Isatin  
**MF:** C<sub>15</sub>H<sub>9</sub>Cl<sub>2</sub>NO<sub>2</sub>  
**FW:** 306.1  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 245, 296 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Apoptosis activator 2 is supplied as a crystalline solid. A stock solution may be made by dissolving the apoptosis activator 2 in the solvent of choice. Apoptosis activator 2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of apoptosis activator 2 in ethanol is approximately 1 mg/ml and 25 mg/ml in DMSO and DMF.

Apoptosis activator 2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, apoptosis activator 2 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Apoptosis activator 2 has a solubility of 0.25 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Apoptosis activator 2 promotes the cytochrome c-dependent oligomerization of Apaf-1 into the mature apoptosome, thus increasing procaspase-9 processing and subsequent caspase-3 activation.<sup>1</sup> This compound has been shown to induce apoptosis in leukemia cells (IC<sub>50</sub> = 4-9 μM) with weak or no effect on normal cell lines or those lacking Apaf-1, caspase-9, or caspase-3 activity (IC<sub>50</sub>s > 40 μM).<sup>1</sup>

### Reference

1. Nguyen, J. T. and Wells, J. A. Direct activation of the apoptosis machinery as a mechanism to target cancer cells. *PNAS* **100**(3), 7533-7538 (2003).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/09/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM