

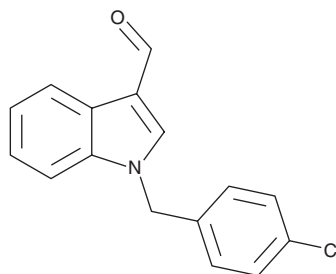
# PRODUCT INFORMATION



## Oncrasin-1

Item No. 25808

CAS Registry No.: 75629-57-1  
Formal Name: 1-[(4-chlorophenyl)methyl]-1H-indole-3-carboxaldehyde  
MF:  $C_{16}H_{12}ClNO$   
FW: 269.7  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{max}$ : 213, 245, 300 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Oncrasin-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the oncrasin-1 in the solvent of choice, which should be purged with an inert gas. Oncrasin-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of oncrasin-1 in these solvents is approximately 5, 17, and 10 mg/ml, respectively.

Oncrasin-1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, oncrasin-1 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Oncrasin-1 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Oncrasin-1 is an anticancer agent that has activity against lung cancer cells expressing constitutively active K-Ras ( $IC_{50}s = 0.25, 1, 1.58, \text{ and } 0.11$  for H460, H2122, A549, and H2887 cells, respectively).<sup>1</sup> It induces nuclear aggregation of atypical protein kinase C (aPKC) PKC $\iota$  and apoptosis in T29Kt1 and H460 cells. Oncrasin-1 (100 mg/kg per day) reduces tumor volume and increases survival time in an H460 mouse xenograft model.

### Reference

1. Guo, W., Wu, S., Liu, J., *et al.* Identification of a small molecule with synthetic lethality for K-Ras and protein kinase C  $\iota$  (PKC $\iota$ ). *Cancer Res.* **68**(18), 7403-7408 (2008).

#### 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, 11/10/2022

#### 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