

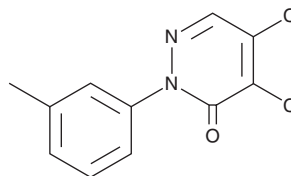
# PRODUCT INFORMATION



## LCS-1

Item No. 35231

**CAS Registry No.:** 41931-13-9  
**Formal Name:** 4,5-dichloro-2-(3-methylphenyl)-3(2H)-pyridazinone  
**Synonym:** Lung Cancer Screen 1  
**MF:** C<sub>11</sub>H<sub>8</sub>Cl<sub>2</sub>N<sub>2</sub>O  
**FW:** 255.1  
**Purity:** ≥98%  
**Supplied as:** A 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

LCS-1 is supplied as a solid. A stock solution may be made by dissolving the LCS-1 in the solvent of choice, which should be purged with an inert gas. LCS-1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LCS-1 in these solvents is approximately 3 and 2 mg/ml, respectively.

### Description

LCS-1 is an inhibitor of superoxide dismutase 1 (SOD1; IC<sub>50</sub> = 1.07 μM).<sup>1</sup> It inhibits the proliferation of NCI H358 non-small cell lung cancer (NSCLC) cells, an effect that can be decreased by overexpression of SOD1 (IC<sub>50</sub> = 0.8 and 3.5 μM, respectively).

### Reference

1. Somwar, R., Erdjument-Bromage, H., Larsson, E., *et al.* Superoxide dismutase 1 (SOD1) is a target for a small molecule identified in a screen for inhibitors of the growth of lung adenocarcinoma cell lines. *Proc. Natl. Acad. Sci. USA* **108**(39), 16375-16380 (2011).

#### 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, 10/06/2022

#### 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](http://WWW.CAYMANCHEM.COM)