

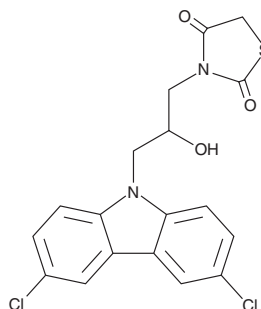
# PRODUCT INFORMATION



**10074-A4**

Item No. 37486

**CAS Registry No.:** 312631-87-1  
**Formal Name:** 3-[3-(3,6-dichloro-9H-carbazol-9-yl)-2-hydroxypropyl]-2,4-thiazolidinedione  
**MF:** C<sub>18</sub>H<sub>14</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>3</sub>S  
**FW:** 409.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 233, 240 nm  
**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

10074-A4 is supplied as a solid. A stock solution may be made by dissolving the 10074-A4 in the solvent of choice, which should be purged with an inert gas. 10074-A4 is soluble in DMSO.

## Description

10074-A4 is an inhibitor of the protein-protein interaction between c-Myc and the c-Myc associated protein Max.<sup>1</sup> It inhibits c-Myc and Max heterodimerization when used at a concentration of 5 μM and reduces the proliferation of Rat1 and TGR-1 rat fibroblasts but not TGR-1 Myc knockout fibroblasts at 32 μM. 10074-A4 (32 μM) reduces the expression of the gene encoding ornithine decarboxylase (ODC), a c-Myc and Max target gene, and induces cell cycle arrest at the G<sub>0</sub>/G<sub>1</sub> phase in Myc-overexpressing Rat1 cells. Preincubation of 10074-A4 (32 μM) with Myc-overexpressing Rat1 cells *in vitro* reduces tumor growth of the same cells in a mouse xenograft model.

## Reference

1. Yin, X., Giap, C., Lazo, J.S., *et al.* Low molecular weight inhibitors of Myc-Max interaction and function. *Oncogene* **22**(40), 6151-6159 (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, 03/28/2023

## 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)