

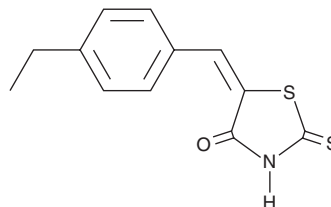
# PRODUCT INFORMATION



**10058-F4**

Item No. 15929

**CAS Registry No.:** 403811-55-2  
**Formal Name:** 5-[(4-ethylphenyl)methylene]-2-thioxo-4-thiazolidinone  
**MF:** C<sub>12</sub>H<sub>11</sub>NOS<sub>2</sub>  
**FW:** 249.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 233, 277, 379 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

10058-F4 is supplied as a crystalline solid. A stock solution may be made by dissolving the 10058-F4 in the solvent of choice, which should be purged with an inert gas. 10058-F4 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 10058-F4 in these solvents is approximately 0.3, 25, and 30, respectively.

10058-F4 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 10058-F4 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 10058-F4 has a solubility of approximately 0.5 in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

The Myc proto-oncogenes interact with Max to form a dimer that regulates gene transcription. The protein c-Myc, in particular, promotes gene expression relevant to cell growth and thus drives cancer. 10058-F4 is a cell-permeable thiazolidinone that inhibits the dimerization of c-Myc and Max at 64 μM, preventing c-Myc-dependent gene expression and cell proliferation.<sup>1</sup> It induces cell cycle arrest, apoptosis, and myeloid differentiation at 100 μM in human acute myeloid leukemia cells.<sup>2</sup> 10058-F4 is rapidly metabolized in mice when given intravenously, limiting its effects on tumors *in vivo*.<sup>3</sup> In addition to c-Myc, 10058-F4 inhibits the nuclear Myc protein, N-Myc, at 50 μM, inducing arrest, apoptosis, and differentiation in neuroblastoma cells overexpressing the gene for N-Myc.<sup>4</sup> This compound can be used to delineate novel actions of Myc proteins, especially those related to lipid and glucose metabolism.<sup>4,5</sup>

## References

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).
2. Huang, M.-J., Cheng, Y.-C., Liu, C.-R., *et al.* A small-molecule c-Myc inhibitor, 10058-F4, induces cell-cycle arrest, apoptosis, and myeloid differentiation of human acute myeloid leukemia. *Exp. Hematol.* **34(11)**, 1480-1489 (2006).
3. Guo, J., Parise, R.A., Joseph, E., *et al.* Efficacy, pharmacokinetics, tissue distribution, and metabolism of the Myc-Max disruptor, 10058-F4 [Z,E]-5-[4-ethylbenzylidene]-2-thioxothiazolidin-4-one, in mice. *Cancer Chemother. Pharmacol.* **63(4)**, 615-625 (2009).
4. Zirath, H., Frenzel, A., Oliynyk, G., *et al.* MYC inhibition induces metabolic changes leading to accumulation of lipid droplets in tumor cells. *Proc. Natl. Acad. Sci. USA* **110(25)**, 10258-10263 (2013).
5. Zhang, P., Metukuri, M.R., Bindom, S.M., *et al.* c-Myc is required for the ChREBP-dependent activation of glucose-responsive genes. *Mol. Endocrinol.* **24(6)**, 1274-1286 (2010).

### 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