

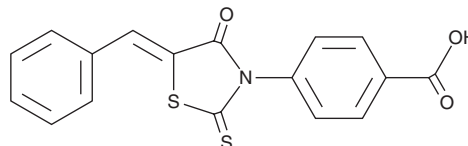
# PRODUCT INFORMATION



## BML-260

Item No. 42442

CAS Registry No.: 101439-76-3  
Formal Name: 4-[4-oxo-5-(phenylmethylene)-2-thioxo-3-thiazolidinyl]-benzoic acid  
MF:  $C_{17}H_{11}NO_3S_2$   
FW: 341.4  
Purity:  $\geq 98\%$   
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

BML-260 is supplied as a crystalline solid. A stock solution may be made by dissolving the BML-260 in the solvent of choice, which should be purged with an inert gas. BML-260 is soluble ( $\geq 10$  mg/ml) in DMSO.

### Description

BML-260 is an inhibitor of dual-specificity phosphatase 22 (DUSP22).<sup>1</sup> *In vivo*, BML-260 (1 mg/kg) increases expression of *Ucp1*, the gene encoding the thermogenic protein uncoupling protein 1 (Ucp1), in white adipocytes and induces thermogenesis in mice.<sup>1</sup> Xenotransplantation of BML-260-expanded hematopoietic stem cells (HSCs) increases immune cell content in immunocompromised mice.<sup>2</sup>

### References

1. Feng, Z., Wei, Y., Zhang, Y., *et al.* Identification of a rhodanine derivative BML-260 as a potent stimulator of UCP1 expression. *Theranostics* **9**(12), 3501-3514 (2019).
2. Albayrak, E., Akgol, S., Turan, R.D., *et al.* BML-260 promotes the growth of cord blood and mobilized peripheral blood hematopoietic stem and progenitor cells with improved reconstitution ability. *J. Cell Biochem.* **123**(12), 2009-2029 (2022).

#### 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, 01/27/2025

#### 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