

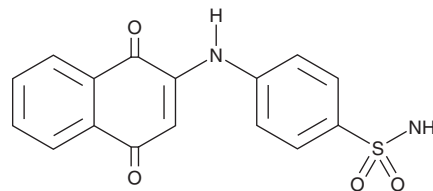
# PRODUCT INFORMATION



## ML-329

Item No. 22481

**CAS Registry No.:** 19992-50-8  
**Formal Name:** 4-[(1,4-dihydro-1,4-dioxo-2-naphthalenyl)amino]-benzenesulfonamide  
**MF:** C<sub>16</sub>H<sub>12</sub>N<sub>2</sub>O<sub>4</sub>S  
**FW:** 328.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 202, 281, 455 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

ML-329 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-329 in the solvent of choice, which should be purged with an inert gas. ML-329 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ML-329 in ethanol is approximately 0.1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

ML-329 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ML-329 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ML-329 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

ML-329 is an inhibitor of the microphthalmia-associated transcription factor (MITF) pathway with an IC<sub>50</sub> value of 1.2 μM for promoter activity of the MITF target gene TRPM-1.<sup>1</sup> It is selectively cytotoxic to the SK-MEL-5 and MALME-3M MITF-dependent melanoma cell lines over A375 MITF-independent cells (IC<sub>50</sub>s = 0.1, 0.7, and 70 μM, respectively).

### Reference

1. Faloon, P.W., Bennion, M., Weiner, W.S., *et al.* A small molecule inhibitor of the MITF molecular pathway. *Probe Reports from the NIH Molecular Libraries Program* (2012).

#### 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, 12/19/2022

#### 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