PRODUCT INFORMATION



ML-334

Item No. 31439

CAS Registry No.: Formal Name:	1432500-66-7 (1S,2R)-2-[[(1S)-1-[(1,3-dihydro- 1,3-dioxo-2H-isoindol-2-yl) methyl]-3,4-dihydro-2(1H)- isoquinolinyl]carbonyl]- cyclohexanecarboxylic acid	
Synonym:	LH601	0 0
MF:	C ₂₆ H ₂₆ N ₂ O ₅	\sim \downarrow \downarrow \downarrow
FW:	446.5	
Purity:	≥98%	
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥2 years	

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

ML-334 is supplied as a solid. A stock solution may be made by dissolving the ML-334 in the solvent of choice, which should be purged with an inert gas. ML-334 is soluble in the organic solvent DMSO at a concentration of approximately 40 mg/ml.

Description

ML-334 is an inhibitor of the protein-protein interaction between the transcription factor Nrf2 and its inhibitor Keap1.¹ It binds to the Keap1 Kelch domain (IC₅₀ = 1 μ M) to induce dissociation and nuclear translocation of Nrf2 and activation of the antioxidant response element (ARE) in a cell-based assay (EC₅₀ = 12 μM).

Reference

1. Hu, L., Magesh, S., Chen, L., et al. Discovery of a small-molecule inhibitor and cellular probe of Keap1-Nrf2 protein-protein interaction. Bioorg. Med. Chem. Lett. 23(10), 3039-3043 (2013).

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.

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, 07/27/2020

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