PRODUCT INFORMATION



KL1333

Item No. 39670

CAS Registry No.:	1800405-30-4	
Formal Name:	2-(1-methylethyl)-3H-	H, I
	naphth[1,2-d]imidazole-4,5-dione	
MF:	$C_{14}H_{12}N_{2}O_{2}$	
FW:	240.3	
Purity:	≥98%	
Supplied as:	A solid	\rightarrow
Storage:	-20°C	ó' `o
Stability:	≥4 years	

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

Laboratory Procedures

KL1333 is supplied as a solid. A stock solution may be made by dissolving the KL1333 in the solvent of choice, which should be purged with an inert gas. KL1333 is soluble (≥10 mg/ml) in DMSO.

Description

KL1333 is a substrate for NAD(P)H:quinone acceptor oxioreductase 1 (NQO1) and a derivative of the quinone β -lapachone (Item No. 15021).¹ It is converted by NQO1 to NAD⁺ and increases the ratio of NAD⁺ to NADH in a cell-free assay, as well as C2C12 and L6 myoblasts, when used at concentrations of 1, 1, or 2 μ M, respectively. KL1333 (1 μ M) increases intracellular ATP levels, as well as decreases intracellular lactate and reactive oxygen species (ROS) levels, in primary fibroblasts isolated from patients with the inborn error of metabolism mitochondrial encephalomyopathy, lactic acidosis, and stroke-like episodes (MELAS). It increases the levels of mitochondrial complex I, also known as NADH dehydrogenase, and mitochondrial complex III, also known as cytochrome bc1 complex, as well as increases the oxygen consumption rate (OCR), in primary fibroblasts isolated from patients with MELAS when used at a concentration of 1 μ M. KL1333 (1 μ M) inhibits apoptosis, the production of mitochondrial ROS, and mitochondrial membrane depolarization induced by the DNA-crosslinking agent cisplatin (Item No. 13119) in primary mouse cochlear organ of Corti cells.²

References

- 1. Seo, K.-S., Kim, J.-H., Min, K.-N., et al. KL1333, a novel NAD⁺ Modulator, improves energy metabolism and mitochondrial dysfunction in MELAS fibroblasts. Front. Neurol. 9, 552 (2018).
- 2. Lee, H.-S., Kim, Y.-R., Lee, I.-K., et al. KL1333, a derivative of β-lapachone, protects against cisplatininduced ototoxicity in mouse cochlear cultures. Biomed. Pharmacother. 126, 110068 (2020).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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, 06/03/2024

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