

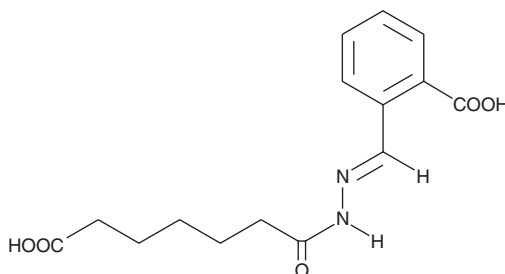
PRODUCT INFORMATION



IDE1

Item No. 13816

CAS Registry No.: 1160927-48-9
Formal Name: 1-[2-[(2-carboxyphenyl)methylene]hydrazide]-heptanedioic acid
MF: C₁₅H₁₈N₂O₅
FW: 306.3
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 285 nm



Laboratory Procedures

For long term storage, we suggest that IDE1 be stored as supplied at -20°C. It should be stable for at least two years.

IDE1 is supplied as a crystalline solid. A stock solution may be made by dissolving the IDE1 in the solvent of choice. IDE1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of IDE1 in these solvents is approximately 0.1, 25, and 30 mg/ml, respectively.

IDE1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, IDE1 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. IDE1 has a solubility of approximately 0.5 mg/ml 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

IDE1 is a small molecule capable of inducing definitive endoderm from embryonic stem cells. It has been shown to induce the differentiation of Sox17+/FoxA2+-expressing pancreatic progenitors from human and mouse embryonic stem cells (EC₅₀ = 125.5 nM *in vitro*) by activating the TGF-β signaling pathway.¹ IDE1-derived endodermal cells injected into E8.75 mouse embryos *ex vivo* have been shown to incorporate into the developing gut tube, contributing to its formation.¹ Furthermore, when treated with either indolactam V (Item No. 14647) or a standard regimen of the growth factor FGF-10, retinoic acid, and hedgehog inhibitors, IDE1-induced endodermal cells can form Pdx1-expressing pancreatic progenitors.¹

Reference

1. Borowiak, M., Maehr, R., Chen, S., *et al.* Small molecules efficiently direct endodermal differentiation of mouse and human embryonic stem cells. *Cell Stem Cell* 4, 348-358 (2009).

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, 09/02/2015

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