

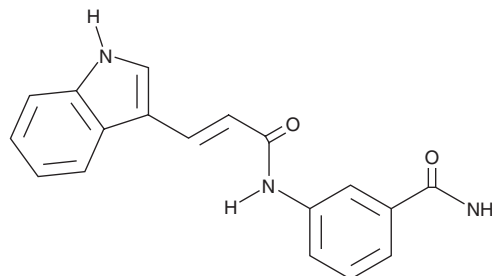
PRODUCT INFORMATION



RSC-133

Item No. 9001839

CAS Registry No.: 1418131-46-0
Formal Name: 3-[[[(2E)-3-(1H-indol-3-yl)-1-oxo-2-propen-1-yl]amino]-benzamide
MF: C₁₈H₁₅N₃O₂
FW: 305.3
Purity: ≥95%
UV/Vis.: λ_{max}: 224, 277, 340, 351 nm
Supplied as: A crystalline 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

RSC-133 is supplied as a crystalline solid. A stock solution may be made by dissolving the RSC-133 in the solvent of choice, which should be purged with an inert gas. RSC-133 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of RSC-133 in these solvents is approximately 5, 20, and 25 mg/ml, respectively.

RSC-133 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RSC-133 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. RSC-133 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

RSC-133 is an indole compound that promotes the reprogramming of human somatic cells to pluripotent stem cells.¹ Specifically, when used at 10 μM with four standard reprogramming factors, it increases the number of human foreskin fibroblasts that express the stem cell marker alkaline phosphatase.¹ RSC-133 induces pluripotency by both down-regulating inducers of cellular senescence and inhibiting DNA methyltransferase and histone deacetylase activities.¹

Reference

1. Lee, J., Xia, Y., Son, M.Y., *et al.* A novel small molecule facilitates the reprogramming of human somatic cells into a pluripotent state and supports the maintenance of an undifferentiated state of human pluripotent stem cells. *Angew. Chem. Int. Ed.* **51(50)**, 12509-12513 (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, 06/26/2020

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