

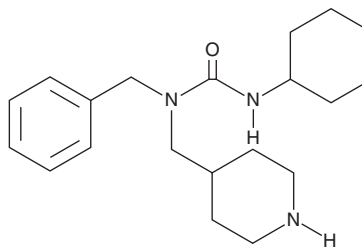
PRODUCT INFORMATION



SRI-011381

Item No. 28914

CAS Registry No.: 1629138-41-5
Formal Name: N'-cyclohexyl-N-(phenylmethyl)-
N-(4-piperidinylmethyl)-urea
MF: C₂₀H₃₁N₃O
FW: 329.5
Purity: ≥98%
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

SRI-011381 is supplied as a crystalline solid. A stock solution may be made by dissolving the SRI-011381 in the solvent of choice, which should be purged with an inert gas. SRI-011381 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SRI-011381 in these solvents is approximately 25, 20, and 15 mg/ml, respectively.

SRI-011381 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SRI-011381 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SRI-011381 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SRI-011381 is an activator of TGF- β signaling.¹ It reduces increases in cell death and the number of dystrophic neurites induced by amyloid- β (1-42) (A β 42; Item No. 20574) in primary mouse embryonic forebrain neurons when used at a concentration of 3 μ M. SRI-011381 (2 and 5 μ M) increases phagocytosis of A β 42 by greater than 20% in J774A.1 and THP-1 macrophages. It increases contextual fear conditioning freezing time and spontaneous alternations in the Y-maze, indicating prevention of memory deficits, in the APP751^{Lon,Swe} transgenic mouse model of Alzheimer's disease when administered at a dose of 10 mg/kg for 10 weeks.

Reference

1. Tanga, M.J. Benzyl urea derivatives for activating TGF- β signaling. *The Board of Trustees of the Leland Stanford Junior University*. **US 2016/0039756 A1** (2016).

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, 12/08/2022

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