

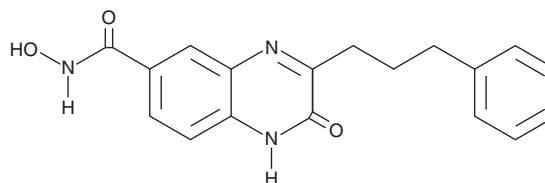
PRODUCT INFORMATION



TCS 2210

Item No. 16142

CAS Registry No.: 1201916-31-5
Formal Name: 1,2-dihydro-N-hydroxy-2-oxo-3-(3-phenylpropyl)-6-quinoxalinecarboxamide
MF: C₁₈H₁₇N₃O₃
FW: 323.3
Purity: ≥95%
Supplied as: A 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

TCS 2210 is supplied as a solid. A stock solution may be made by dissolving the TCS 2210 in the solvent of choice. TCS 2210 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 5 mg/ml.

Description

TCS 2210 is a small molecule inducer of neuronal differentiation.¹ It increases expression of the neuronal markers β -III tubulin and neuron-specific enolase (NSE) and induces neurite outgrowth in a population of PC12 neuronal precursor-like cells. TCS 2210 (20 μ M per day for two days) converts >95% of a rat mesenchymal stem cell (MSC) population to a neuronal phenotype that exhibits potassium outward currents and increased expression of β -III tubulin and NSE as well as the cholinergic genes *CHRNA2*, *CHRNAB*, and *CHRM4*.

Reference

1. Kim, N.R., Kang, S.K., Ahn, H.H., et al. Discovery of a new and efficient small molecule for neuronal differentiation from mesenchymal stem cell. *J. Med. Chem.* **52(24)**, 7931-7933 (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, 12/05/2022

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