

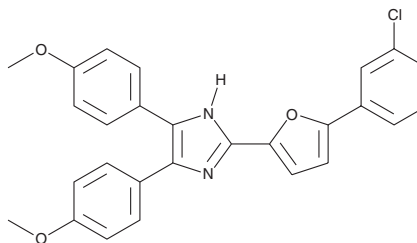
PRODUCT INFORMATION



Neurodazine

Item No. 13224

CAS Registry No.: 937807-66-4
Formal Name: 2-[5-(3-chlorophenyl)-2-furanyl]-4,5-bis(4-methoxyphenyl)-1H-imidazole
MF: C₂₇H₂₁ClN₂O₃
FW: 456.9
Purity: ≥98%
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

Neurodazine is supplied as a crystalline solid. A stock solution may be made by dissolving the neurodazine in the solvent of choice, which should be purged with an inert gas. Neurodazine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of neurodazine in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

Neurodazine induces neuronal differentiation in skeletal muscle cells.¹ In mouse C2C12 myoblasts, neurodazine (2 μM for 7 days) upregulated the expression of neuron-specific enolase (NSE), neurofilament 200 (NF200), synapsin, and other neuronal markers. Neurodazine similarly upregulated NSE and NF200 in human skeletal muscle fibers, indicating efficacy in mature muscle as well as in myoblasts.

Reference

1. Williams, D.R., Lee, M.-R., Song, Y.-A., *et al.* Synthetic small molecules that induce neurogenesis in skeletal muscle. *J. Am. Chem. Soc.* **129**, 9258-9259 (2007).

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, 06/25/2020

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