

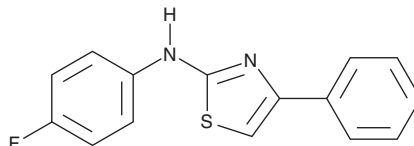
PRODUCT INFORMATION



UCLA gp130 2

Item No. 35551

CAS Registry No.: 339303-87-6
Formal Name: N-(4-fluorophenyl)-4-phenyl-2-thiazolamine
MF: $C_{15}H_{11}FN_2S$
FW: 270.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 243, 278 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

UCLA gp130 2 is supplied as a solid. A stock solution may be made by dissolving the UCLA gp130 2 in the solvent of choice, which should be purged with an inert gas. UCLA gp130 2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of UCLA gp130 2 in these solvents is approximately 5, 3, and 1 mg/ml, respectively.

Description

UCLA gp130 2 is a humanin peptide mimetic.¹ It inhibits NMDA-induced lactate dehydrogenase (LDH) release, a marker of cytotoxicity, in primary mouse hippocampal neurons when used at a concentration of 10 μM , an effect that can be reversed by the gp130 inhibitor SC144 (Item No. 29714).

Reference

1. Alam, M.P., Bilousova, T., Spilman, P., *et al.* A small molecule mimetic of the humanin peptide as a candidate for modulating NMDA-Induced neurotoxicity. *ACS Chem. Neurosci.* **9**(3), 462-468 (2018).

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, 12/12/2022

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