

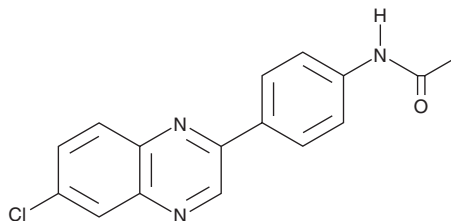
PRODUCT INFORMATION



CA77.1

Item No. 37300

CAS Registry No.: 2412270-22-3
Formal Name: N-[4-(6-chloro-2-quinoxalinyloxy)]-acetamide
MF: C₁₆H₁₂ClN₃O
FW: 297.7
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 269, 302, 363 nm
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

CA77.1 is supplied as a solid. A stock solution may be made by dissolving the CA77.1 in the solvent of choice, which should be purged with an inert gas. CA77.1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CA77.1 in these solvents is approximately 1 and 2 mg/ml, respectively.

Description

CA77.1 is an inducer of chaperone-mediated autophagy (CMA).¹ It induces the formation of autophagic lysosomes in NIH3T3 cells expressing a CMA reporter when used at concentrations of 10, 20, and 30 μM. CA77.1 (20 μM) induces autophagic flux in NIH3T3 cells. *In vivo*, CA77.1 (30 mg/kg) reduces locomotor hyperactivity to control levels and the number of neurons containing pathogenic tau conformations in the hippocampus, amygdala, and piriform cortex in a PS19 transgenic mouse model of Alzheimer's disease.

Reference

1. Bourdenx, M., Martin-Segura, A., Scrivo, A., *et al.* Chaperone-mediated autophagy prevents collapse of the neuronal metastable proteome. *Cell* **184**(10), 2696-2714.e25 (2021).

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, 02/10/2023

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