

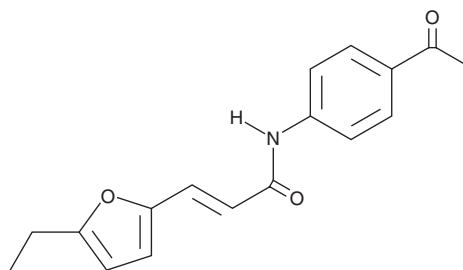
PRODUCT INFORMATION



Heclin

Item No. 26139

CAS Registry No.: 890605-54-6
Formal Name: N-(4-acetylphenyl)-3-(5-ethyl-2-furanyl)-2-propenamide
MF: C₁₇H₁₇NO₃
FW: 283.3
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 342 nm
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

Heclin is supplied as a crystalline solid. A stock solution may be made by dissolving the heclin in the solvent of choice, which should be purged with an inert gas. Heclin is soluble in organic solvents such as ethanol and DMSO. The solubility of heclin in these solvents is approximately 10 and 20 mg/ml, respectively.

Description

Heclin is an inhibitor of HECT E3 ubiquitin ligases (IC₅₀s = 6.8, 6.3, and 6.9 μM for Smurf2, Nedd4, and WWP1 HECT ligase domains, respectively).¹ It inhibits autoubiquitination of Smurf2 in HEK293 cells expressing the human enzyme (IC₅₀ = 9 μM) and reduces ubiquitination of Dishevelled 2 in HEK293 cells expressing Smurf2, Nedd4, or WWP2. Heclin is cytotoxic to HEK293 cells (IC₅₀ = 45 μM).

Reference

1. Mund, T., Lewis, M.J., Maslen, S., *et al.* Peptide and small molecule inhibitors of HECT-type ubiquitin ligases. *Proc. Natl. Acad. Sci. U.S.A.* **111**(47), 16736-16741 (2014).

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, 12/02/2022

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