

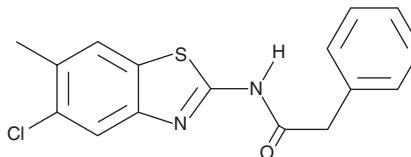
PRODUCT INFORMATION



LH 846

Item No. 17686

CAS Registry No.: 639052-78-1
Formal Name: N-(5-chloro-6-methyl-2-benzothiazolyl)-benzeneacetamide
MF: C₁₆H₁₃ClN₂OS
FW: 316.8
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 233, 274, 296, 307 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

LH 846 is supplied as a crystalline solid. A stock solution may be made by dissolving the LH 846 in the solvent of choice, which should be purged with an inert gas. LH 846 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LH 846 in these solvents is approximately 5 mg/ml.

Description

LH 846 is a benzothiazole analog that selectively inhibits casein kinase (CK) 1δ (IC₅₀ = 290 nM).¹ It inhibits CK1ε and CK1α at much higher concentrations (IC₅₀s = 1.3 and 2.5 μM, respectively) but does not inhibit CK2.¹ LH 846 has been shown to inhibit CK1δ-dependent phosphorylation and degradation of PER1 protein and to lengthen the circadian period in U2OS cells.¹

Reference

1. Lee, J.W., Hirota, T., Peters, E.C., M., *et al.* A small molecule modulates circadian rhythms through phosphorylation of the period protein. *Angew Chem. Int. Ed. Engl.* **50(45)**, 10608-10611 (2011).

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, 11/08/2022

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