

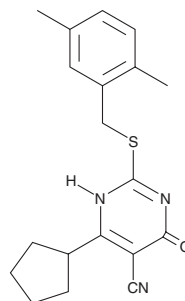
# PRODUCT INFORMATION



**HJC0197**

Item No. 19092

**CAS Registry No.:** 1383539-73-8  
**Formal Name:** 4-cyclopentyl-2-[[[(2,5-dimethylphenyl)methyl]thio]-1,6-dihydro-6-oxo-5-pyrimidinecarbonitrile  
**MF:** C<sub>19</sub>H<sub>21</sub>N<sub>3</sub>OS  
**FW:** 339.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 203, 312 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

HJC0197 is supplied as a crystalline solid. A stock solution may be made by dissolving the HJC0197 in the solvent of choice. HJC0197 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of HJC0197 in these solvents is approximately 5 and 10 mg/ml, respectively.

HJC0197 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HJC0197 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. HJC0197 has a solubility of approximately 0.2 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

Exchange protein activated by cAMP (Epac) proteins mediate cAMP signaling independent of protein kinase A (PKA). HJC0197 is a cell-permeable inhibitor of Epac1 and Epac2 (IC<sub>50</sub> = 5.9 μM for Epac2).<sup>1</sup> It inhibits Epac1-mediated Rap1-GDP exchange activity at 25 μM, but has no effect on cAMP-induced type I and II PKA activity at this concentration.<sup>1</sup> Pretreatment of HEK293 cells expressing either Epac1 or Epac2 with 10 μM HJC0197 completely blocks Epac-mediated phosphorylation of Akt.<sup>1</sup> HJC0197 has been used to study the role of Epac signaling in chondrogenesis in chicken micromass cultures.<sup>2</sup>

## References

1. Chen, H., Tsalkova, T., Mei, F.C., *et al.* 5-Cyano-6-oxo-1,6-dihydro-pyrimidines as potent antagonists targeting exchange proteins directly activated by cAMP. *Bioorg. Med. Chem. Lett.* **22**(12), 4038-4043 (2012).
2. Juhász, T., Matta, C., Somogyi, C., *et al.* Mechanical loading stimulates chondrogenesis via the PKA/CREB-Sox9 and PP2A pathways in chicken micromass cultures. *Cell Signal.* **26**, 468-482 (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, 12/19/2022

## 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](http://WWW.CAYMANCHEM.COM)