

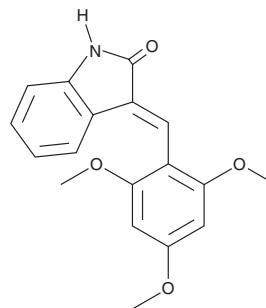
# PRODUCT INFORMATION



## IC261

Item No. 17260

**CAS Registry No.:** 186611-52-9  
**Formal Name:** 1,3-dihydro-3-[(2,4,6-trimethoxyphenyl)methylene]-2H-indol-2-one  
**Synonym:** SU5607  
**MF:** C<sub>18</sub>H<sub>17</sub>NO<sub>4</sub>  
**FW:** 311.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 255, 358 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

IC261 is supplied as a crystalline solid. A stock solution may be made by dissolving the IC261 in the solvent of choice. IC261 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of IC261 in ethanol is approximately 0.5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

IC261 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, IC261 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. IC261 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

IC261 is a reversible, ATP-competitive inhibitor of casein kinase 1 (CK1) that inhibits CK1δ and CK1ε (IC<sub>50</sub> = ~1 μM for both), as well as CK1α (IC<sub>50</sub> = 16 μM).<sup>1</sup> It is at least 100-fold less effective against PKA, p34<sup>cdc2</sup>, and p55<sup>fyn</sup>.<sup>1</sup> IC261, at 1 μM, inhibits cytokinesis in primary mouse embryo fibroblasts.<sup>2</sup> IC261 is used to elucidate the role of CK1 in cells and in whole organisms.<sup>3-5</sup>

### References

1. Mashhoon, N., DeMaggio, A.J., Tereshko, V., *et al.* Crystal structure of a conformation-selective casein kinase-1 inhibitor. *J. Biol. Chem.* **275**(26), 20052-20060 (2000).
2. Behrend, L., Milne, D.M., Stöter, M., *et al.* IC261, a specific inhibitor of the protein kinases casein kinase 1-delta and -epsilon, triggers the mitotic checkpoint and induces p53-dependent postmitotic effects. *Oncogene* **19**(47), 5303-5313 (2000).
3. van Ooijen, G., Martin, S.F., Barrios-Llerena, M.E., *et al.* Functional analysis of the rodent CK1<sup>tau</sup> mutation in the circadian clock of a marine unicellular alga. *BMC Cell Biol.* **14**, (2013).
4. Rachidi, N., Taly, J.F., Durieu, E., *et al.* Pharmacological assessment defines *Leishmania donovani* casein kinase 1 as a drug target and reveals important functions in parasite viability and intracellular infection. *Antimicrob. Agents Chemother.* **58**(3), 1501-1515 (2014).
5. Kurihara, T., Sakurai, E., Toyomoto, M., *et al.* Alleviation of behavioral hypersensitivity in mouse models of inflammatory pain with two structurally different casein kinase 1 (CK1) inhibitors. *Mol. Pain* **10**, (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, 10/13/2022

#### 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