

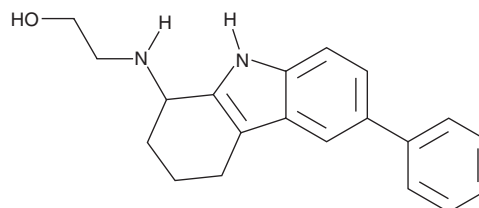
# PRODUCT INFORMATION



## CASIN

Item No. 17694

**CAS Registry No.:** 425399-05-9  
**Formal Name:** 2-[(2,3,4,9-tetrahydro-6-phenyl-1H-carbazol-1-yl)amino]-ethanol  
**Synonym:** Pirl1-related Compound 2  
**MF:** C<sub>20</sub>H<sub>22</sub>N<sub>2</sub>O  
**FW:** 306.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 255 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

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

### Description

CASIN is an indole-based small molecule that inhibits activation of the Cdc42 GTPase (IC<sub>50</sub> = 2 μM for inhibiting actin assembly in PIP<sub>2</sub>-stimulated actin polymerization assays).<sup>1</sup> This compound has been used to explore the role of Cdc42 in the division, survival, and differentiation of hematopoietic stem cells.<sup>2,3</sup>

### References

1. Peterson, J.R., Lebensohn, A.M., Pelish, H.E., *et al.* Biochemical suppression of small molecule inhibitors: A new strategy to identify inhibitor targets and signaling pathway components. *Chem. Biol.* **13(4)**, 443-452 (2006).
2. Sakamori, R., Das, S., Yu, S., *et al.* Cdc42 and Rab8a are critical for intestinal stem cell division, survival, and differentiation in mice. *J. Clin. Invest.* **122(3)**, 1052-1065 (2012).
3. Florian, M.C., Dörr, K., Niebel, A., *et al.* Cdc42 activity regulates hematopoietic stem cell aging and rejuvenation. *Cell Stem Cell* **10(5)**, 520-530 (2012).

#### 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