

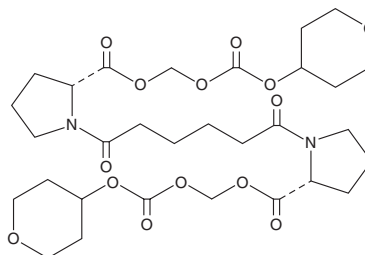
# PRODUCT INFORMATION



**GSK3039294**

Item No. 36559

**CAS Registry No.:** 1819986-22-5  
**Formal Name:** 1,1'-(1,6-dioxo-1,6-hexanediyl)bis-D-proline, 2,2'-bis[[[(tetrahydro-2H-pyran-4-yl)oxy]carbonyl]oxy]methyl] ester  
**MF:** C<sub>30</sub>H<sub>44</sub>N<sub>2</sub>O<sub>14</sub>  
**FW:** 656.7  
**Purity:** ≥95%  
**Supplied as:** A 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

GSK3039294 is supplied as a solid. A stock solution may be made by dissolving the GSK3039294 in the solvent of choice, which should be purged with an inert gas. GSK3039294 is sparingly soluble (1-10 mg/ml) in DMSO.

## Description

GSK3039294 is a prodrug form of CPHPC (Item No. 75500), an inhibitor of the protein-protein interaction between serum amyloid P (SAP) and amyloid fibrils.<sup>1</sup> It is converted to CPHPC by hydrolysis via a monoester intermediate in rat, canine, and human liver microsomes.

## Reference

- Richards, D., Bamford, M., Liefwaard, L., *et al.* Identification, preclinical profile, and clinical proof of concept of an orally bioavailable pro-drug of miridesap. *Br. J. Pharmacol.* **177**(8), 1853-1864 (2020).

### 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, 05/07/2024

## 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)