

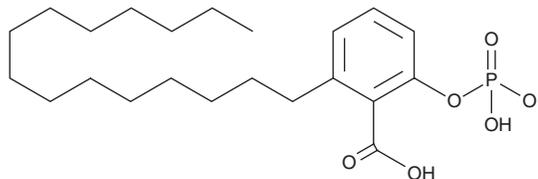
# PRODUCT INFORMATION



## Phospho-Ginkgolic Acid

Item No. 29638

**CAS Registry No.:** 2281761-58-6  
**Formal Name:** 2-pentadecyl-6-(phosphonoxy)-benzoic acid  
**Synonym:** Ginkgolic Acid 2-Phosphate  
**MF:** C<sub>22</sub>H<sub>37</sub>O<sub>6</sub>P  
**FW:** 428.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 244 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Phospho-ginkgolic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the phospho-ginkgolic acid in the solvent of choice, which should be purged with an inert gas. Phospho-ginkgolic acid is soluble in organic solvents such as ethanol and DMSO.

### Description

Phospho-ginkgolic acid is an agonist of sphingosine-1-phosphate receptor 1 (S1P<sub>1</sub>).<sup>1</sup> Phospho-ginkgolic acid (10 μM) induces S1P<sub>1</sub> receptor internalization and increases ERK phosphorylation in a concentration-dependent manner in CHO cells expressing S1P<sub>1</sub>.

### Reference

1. Swamy, M.M.M., Murai, Y., Ohno, Y., *et al.* Structure-inspired design of a sphingolipid mimic sphingosine-1-phosphate receptor agonist from a naturally occurring sphingomyelin synthase inhibitor. *Chem. Commun.* **54(90)**, 12758-12761 (2018).

**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, 01/13/2020

### 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