

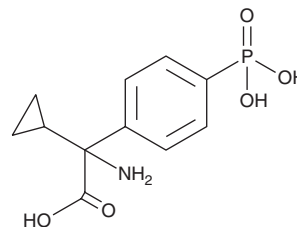
# PRODUCT INFORMATION



## CPPG

Item No. 34362

**CAS Registry No.:** 183364-82-1  
**Formal Name:**  $\alpha$ -amino- $\alpha$ -cyclopropyl-4-phosphono-benzeneacetic acid  
**Synonym:**  $\alpha$ -Cyclopropyl-4-phosphonophenylglycine  
**MF:** C<sub>11</sub>H<sub>14</sub>NO<sub>5</sub>P  
**FW:** 271.2  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

CPPG is supplied as a solid. A stock solution may be made by dissolving the CPPG in the solvent of choice, which should be purged with an inert gas. CPPG is soluble in 1 M sodium hydroxide.

### Description

CPPG is an antagonist of group II and group III metabotropic glutamate receptors (mGluRs).<sup>1</sup> It reverses inhibition of forskolin-induced cAMP accumulation induced by the mGluR group II selective agonist L-CCG-1 or the mGluR group III-selective agonist L-AP4 (Item No. 14538) in rat cerebrocortical slices (IC<sub>50</sub>s = 46.2 and 2.2 nM, respectively). Intrathecal administration of CPPG reduces formalin-induced flinching in a rat model of inflammatory pain.<sup>2</sup>

### References

1. Toms, N.J., Jane, D.E., Kemp, M.C., *et al.* The effects of (RS)- $\alpha$ -cyclopropyl-4-phosphonophenylglycine ((RS)-CPPG), a potent and selective metabotropic glutamate receptor antagonist. *Br. J. Pharmacol.* **119**(5), 851-854 (1996).
2. Okubo, M., Yamanaka, H., Kobayashi, K., *et al.* Differential expression of mGluRs in rat spinal dorsal horns and their modulatory effects on nocifensive behaviors. *Mol. Pain* **15**, 1-13 (2019).

#### 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, 11/22/2022

#### 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