

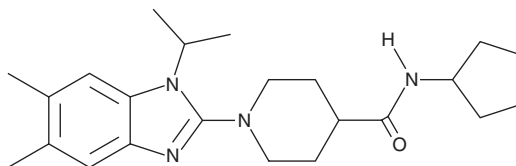
# PRODUCT INFORMATION



**CAY10678**

Item No. 15129

**CAS Registry No.:** 1268709-57-4  
**Formal Name:** N-cyclopentyl-1-[5,6-dimethyl-1-(1-methylethyl)-1H-benzimidazol-2-yl]-4-piperidinecarboxamide  
**Synonym:** mPGES-1 Inhibitor III  
**MF:** C<sub>23</sub>H<sub>34</sub>N<sub>4</sub>O  
**FW:** 382.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 215, 250, 292 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

CAY10678 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10678 in the solvent of choice. CAY10678 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of CAY10678 in these solvents is approximately 25, 5, and 15 mg/ml, respectively.

CAY10678 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10678 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. CAY10678 has a solubility of approximately 0.1 mg/ml in a 1:8 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

Microsomal prostaglandin E synthase-1 (mPGES-1) converts the COX product PGH<sub>2</sub> into the biologically active PGE<sub>2</sub>.<sup>1</sup> Like COX-2, the expression of mPGES-1 is induced in response to pro-inflammatory mediators, including LPS, IL-1β, and TNF-α.<sup>2</sup> CAY10678 is a benzimidazole that potently inhibits human and rat recombinant mPGES-1 (IC<sub>50</sub> = 0.09 and 0.9 μM, respectively).<sup>3</sup> It has minimal effects on COX-1, COX-2, PGIS, and hematopoietic PGDS at 50 μM but reduces lipocalin-type PGDS activity by 60% at this concentration.<sup>3</sup> CAY10678 dose-dependently blocks PGE<sub>2</sub> synthesis in isolated cells and whole blood treated with LPS or IL-1β.<sup>3</sup> It also dose-dependently reduces PGE<sub>2</sub> synthesis and cell recruitment during inflammation in mice.<sup>3</sup>

## References

1. Jakobsson, P.-J., Thorén, S., Morgenstern, R., *et al.* Identification of human prostaglandin E synthase: A microsomal, glutathione-dependent, inducible enzyme, constituting a potential novel drug target. *Proc. Natl. Acad. Sci. USA* **96**, 7220-7225 (1999).
2. Stichtenoth, D.O., Thorén, S., Bian, H., *et al.* Microsomal prostaglandin E synthase is regulated by proinflammatory cytokines and glucocorticoids in primary rheumatoid synovial cells. *J. Immunol.* **167**, 469-474 (2001).
3. Leclerc, P., Idborg, H., Spahiu, L., *et al.* Characterization of a human and murine mPGES-1 inhibitor and comparison to mPGES-1 genetic deletion in mouse models of inflammation. *Prostaglandins Other Lipid Mediat.* (2013).

### 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, 02/01/2024

## 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