

PRODUCT INFORMATION

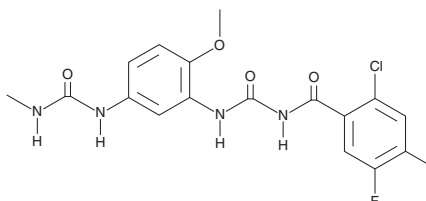


Glycogen Phosphorylase Inhibitor

Item No. 17578

CAS Registry No.: 648926-15-2
Formal Name: 2-chloro-4,5-difluoro-N-[[[2-methoxy-5-[[[(methylamino)carbonyl]amino]phenyl]amino]carbonyl]-benzamide

Synonym: GPI
MF: C₁₇H₁₅ClF₂N₄O₄
FW: 412.8
Purity: ≥98%
UV/Vis.: λ_{max}: 238, 302 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

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

Description

Glycogen phosphorylase in the liver, muscle, and brain initiate glycogenolysis by releasing glucose-1-phosphate from glycogen. Glycogen phosphorylase inhibitor is a cell-permeable acyl urea first identified as an inhibitor of human liver glycogen phosphorylase (IC₅₀ = 53 nM).¹ It blocks glucagon-induced hepatic glycogenolysis *in vivo*.¹ Glycogen phosphorylase inhibitor has been used to study glycogen utilization in human liver HepG2 cells, retinal explants, and human T lymphocyte Kit 225 cells.²⁻⁴

References

1. Klabunde, T., Wendt, K.U., Kadereit, D., *et al.* Acyl ureas as human liver glycogen phosphorylase inhibitors for the treatment of type 2 diabetes. *J. Med. Chem.* **48**(20), 6178-6193 (2005).
2. Cheng, A., Zhang, M., Gentry, M.S., *et al.* A role for AGL ubiquitination in the glycogen storage disorders of Lafora and Cori's disease. *Genes Dev.* **21**(19), 2399-2409 (2007).
3. Agathocleous, M., Love, N.K., Randlett, O., *et al.* Metabolic differentiation in retinal cells. *Nat. Cell Biol.* **14**(8), 859-864 (2012).
4. Arrizabalaga, O., Lacerda, H.M., Zubiaga, A.M., *et al.* Rac1 protein regulates glycogen phosphorylase activation and controls interleukin (IL)-2-dependent T cell proliferation. *J. Biol. Chem.* **287**(15), 11878-11890 (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, 12/12/2022

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