

PRODUCT INFORMATION



CAY10793

Item No. 37436

Formal Name: 7-fluoro-2-(((5-methoxy-1H-benzo[d]imidazol-2-yl)thio)methyl)quinazolin-4(3H)-one

MF: C₁₇H₁₃FN₄O₂S

FW: 356.4

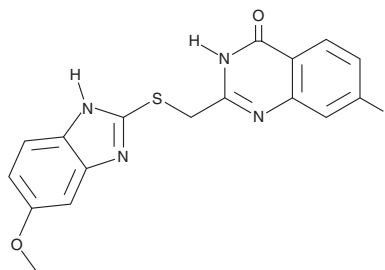
Purity: ≥98%

UV/Vis.: λ_{max}: 298 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

CAY10793 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10793 in the solvent of choice, which should be purged with an inert gas. CAY10793 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CAY10793 in these solvents is approximately 10 mg/ml.

Description

CAY10793 is a quinazolinone-containing compound that is structurally similar to other quinazolinone-containing inhibitors of ENPP1.¹

Reference

1. Wang, X., Lu, X., Yan, D., *et al.* Development of novel ecto-nucleotide pyrophosphatase/phosphodiesterase 1 (ENPP1) inhibitors for tumor immunotherapy. *Int. J. Mol. Sci.* **23(13)**, 7104 (2022).

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, 03/26/2024

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