

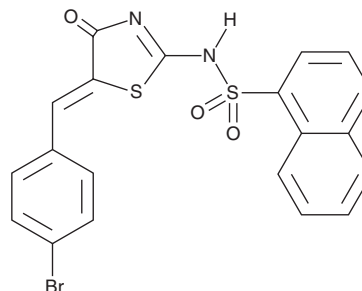
PRODUCT INFORMATION



Pitstop2

Item No. 23885

CAS Registry No.: 1332879-52-3
Formal Name: N-[5-[(4-bromophenyl)methylene]-4,5-dihydro-4-oxo-2-thiazolyl]-1-naphthalenesulfonamide
MF: C₂₀H₁₃BrN₂O₃S₂
FW: 473.4
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 212, 222, 344 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

Pitstop2 is supplied as a crystalline solid. A stock solution may be made by dissolving the pitstop2 in the solvent of choice, which should be purged with an inert gas. Pitstop2 is soluble in the organic solvent DMSO at a concentration of approximately 20 mg/ml.

Description

Pitstop2 is an inhibitor of the interaction between the clathrin terminal domain and amphiphysin (IC₅₀ = 12 μM).¹ It inhibits clathrin-mediated transferrin uptake in HeLa and U2OS cells (IC₅₀s = 12-15 and 9.7 μM, respectively). Pitstop2 inhibits HIV-1 entry and infectivity in HeLa reporter cell lines. It also inhibits clathrin-independent endocytosis of CD44, CD98, and CD147 in HeLa cells.²

References

1. von Kleist, L., Stahlschmidt, W., Bulut, H., *et al.* Role of the clathrin terminal domain in regulating coated pit dynamics revealed by small molecule inhibition. *Cell* **146(3)**, 471-484 (2012).
2. Dutta, D., Williamson, C.D., Cole, N.B., *et al.* Pitstop 2 is a potent inhibitor of clathrin-independent endocytosis. *PLoS One* **7(9)**, e45799 (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, 11/03/2022

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