PRODUCT INFORMATION



WS3

Item No. 17667

| CAS Registry No.: | 1421227-52-2 | |
|-------------------|---|------------------|
| Formal Name: | N-[6-[4-[[[[4-[(4-methyl-1-piperazinyl) | |
| | methyl]-3-(trifluoromethyl)phenyl] | ∠ ^N ∖ |
| | amino]carbonyl]amino]phenoxy]-4- | |
| | pyrimidinyl]-cyclopropanecarboxamide | A N |
| MF: | $C_{28}H_{30}F_3N_7O_3$ | A_{0} |
| FW: | 569.6 | |
| Purity: | ≥95% | |
| UV/Vis.: | λ _{may} : 264 nm | |
| Supplied as: | A crystalline solid | |
| Storage: | -20°C | H |
| Stability: | ≥4 years | |
| | | |

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

Laboratory Procedures

WS3 is supplied as a crystalline solid. A stock solution may be made by dissolving the WS3 in the solvent of choice, which should be purged with an inert gas. WS3 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of WS3 in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

WS3 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, WS3 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. WS3 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

WS3 is a non-specific proliferative molecule that modulates the activity of Erb3 binding protein-1 and the IkB kinase pathway.¹ It has been used to mediate proliferation of primary retinal pigment epithelial cells ex vivo in order to provide a renewable source of cells for transplantation in a model of retinal degeneration.²

References

- 1. Shen, W., Tremblay, M. S., Deshmukh, V. A., et al. Small-molecule inducer of β cell proliferation identified by highthroughput screening. J. Am. Chem. Soc. 135(5), 1669-1672 (2013).
- 2. Swoboda, J.G., Elliott, J., Deshmukh, V., et al. Small molecule mediated proliferation of primary retinal pigment epithelial cells. ACS Chem Biol. 8(7), 1407-1411 (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.

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, 11/10/2022

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