

PRODUCT INFORMATION



WAY-316606

Item No. 26324

CAS Registry No.: 915759-45-4
Formal Name: 5-(phenylsulfonyl)-N-4-piperidinyl-2-(trifluoromethyl)-benzenesulfonamide

MF: C₁₈H₁₉F₃N₂O₄S₂
FW: 448.5

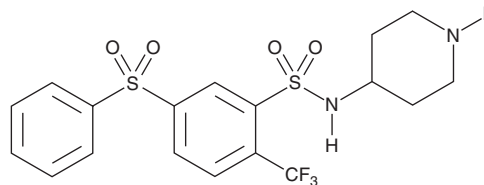
Purity: ≥98%

UV/Vis.: λ_{max}: 239 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

WAY-316606 is supplied as a crystalline solid. A stock solution may be made by dissolving the WAY-316606 in the solvent of choice, which should be purged with an inert gas. WAY-316606 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of WAY-316606 in these solvents is approximately 3 mg/ml.

WAY-316606 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, WAY-316606 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. WAY-316606 has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

WAY-316606 is an inhibitor of secreted frizzled-related protein 1 (sFRP1; IC₅₀ = 0.5 μM).¹ It increases Wnt signaling in U2OS osteosarcoma cells (EC₅₀ = 0.65 μM). WAY-316606 increases total bone area in a murine calvarial organ culture assay when used at concentrations greater than 0.1 nM.²

References

1. Moore, W.J., Kern, J.C., Bhat, R., *et al.* Modulation of Wnt signaling through inhibition of secreted frizzled-related protein I (sFRP-1) with N-substituted piperidinyl diphenylsulfonyl sulfonamides: Part II. *Bioorg. Med. Chem.* **18**(1), 190-201 (2010).
2. Bodine, P.V., Stauffer, B., Ponce-de-Leon, H., *et al.* A small molecule inhibitor of the Wnt antagonist secreted frizzled-related protein-1 stimulates bone formation. *Bone* **44**(6), 1063-1068 (2009).

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, 03/16/2023

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