

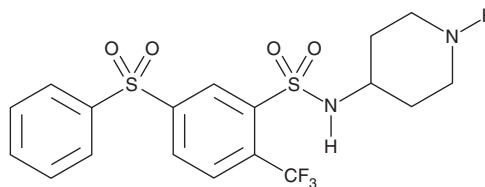
# PRODUCT INFORMATION



**WAY-316606**

Item No. 45425

**CAS Registry No.:** 915759-45-4  
**Formal Name:** 5-(phenylsulfonyl)-N-4-piperidinyl-2-(trifluoromethyl)-benzenesulfonamide  
**MF:** C<sub>18</sub>H<sub>19</sub>F<sub>3</sub>N<sub>2</sub>O<sub>4</sub>S<sub>2</sub>  
**FW:** 448.5  
**Purity:** ≥98%  
**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 sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

## Description

WAY-316606 is an inhibitor of secreted frizzled-related protein 1 (sFRP1; IC<sub>50</sub> = 0.5 μM).<sup>1</sup> It increases Wnt signaling in U2OS osteosarcoma cells (EC<sub>50</sub> = 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.<sup>2</sup>

## 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, 02/23/2026

## 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