

# PRODUCT INFORMATION

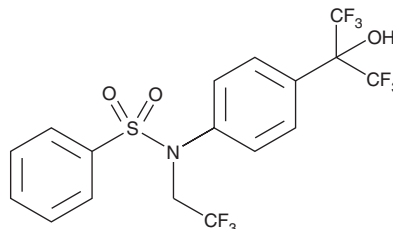


**T0901317**

Item No. 71810

**CAS Registry No.:** 293754-55-9  
**Formal Name:** N-(2,2,2-trifluoroethyl)-N-[4-[2,2,2-trifluoro-1-hydroxy-1-(trifluoromethyl)ethyl]phenyl]-benzenesulfonamide

**MF:** C<sub>17</sub>H<sub>12</sub>F<sub>9</sub>NO<sub>3</sub>S  
**FW:** 481.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 227, 265, 272 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

T0901317 is supplied as a crystalline solid. A stock solution may be made by dissolving the T0901317 in the solvent of choice, which should be purged with an inert gas. T0901317 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of T0901317 in ethanol is approximately 50 mg/ml and approximately 100 mg/ml in DMSO and DMF.

T0901317 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, T0901317 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. T0901317 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

T0901317 is a potent and selective agonist for both LXR $\alpha$  and LXR $\beta$ , with an EC<sub>50</sub> of about 50 nM.<sup>1</sup> T0901317 acting through LXR and in concert with its RXR heterodimerization partner induces the expression of the ABCA1 reverse cholesterol transporter. This acts to increase the efflux of cholesterol from enterocytes and thus inhibit the overall absorption of cholesterol.

## Reference

1. Repa, J.J., Turley, S.D., Lobaccaro, J.-M.A., *et al.* Regulation of absorption and ABC1-mediated efflux of cholesterol by RXR heterodimers. *Science* **289**, 1524-1529 (2000).

### 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, 04/08/2024

## 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