

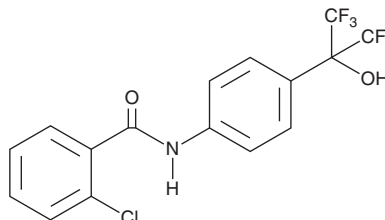
# PRODUCT INFORMATION



**SR 0987**

Item No. 19503

**CAS Registry No.:** 303126-97-8  
**Formal Name:** 2-chloro-N-[4-[2,2,2-trifluoro-1-hydroxy-1-(trifluoromethyl)ethyl]phenyl]-benzamide  
**MF:** C<sub>16</sub>H<sub>10</sub>ClF<sub>6</sub>NO<sub>2</sub>  
**FW:** 397.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 255 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

SR 0987 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 0987 in the solvent of choice, which should be purged with an inert gas. SR 0987 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SR 0987 in these solvents is approximately 30, 15, and 20 mg/ml, respectively.

SR 0987 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SR 0987 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SR 0987 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

SR 0987 is an agonist of the T cell-specific isoform of ROR $\gamma$  (ROR $\gamma$ t, retinoic acid receptor-related orphan receptor- $\gamma$ t), inducing reporter gene expression with an EC<sub>50</sub> value of ~800 nM.<sup>1</sup> Stimulation of mouse EL4 T lymphocytes with SR 0987, following activation with PMA (Item No. 10008014) and ionomycin (Item No. 10004974), increases expression of the ROR $\gamma$ t target IL-17, while expression of the programmed cell death protein PD-1 is decreased and granzyme B is unchanged.<sup>1</sup> Decreased cell surface PD-1 protein, assessed by flow cytometry, is observed in a variety of T cells treated with SR 0987.<sup>1</sup>

## Reference

1. Chang, M.R., Dharmarajan, V., Doebelin, C., *et al.* Synthetic ROR $\gamma$ t agonists enhance protective immunity. *ACS Chem. Biol.* **11**(4), 1012-1018 (2016).

### 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, 12/02/2022

## 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