

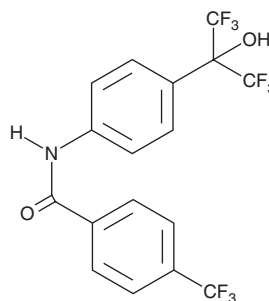
PRODUCT INFORMATION



SR 1078

Item No. 16503

CAS Registry No.: 1246525-60-9
Formal Name: N-[4-[2,2,2-trifluoro-1-hydroxy-1-(trifluoromethyl)ethyl]phenyl]-4-(trifluoromethyl)-benzamide
MF: C₁₇H₁₀F₉NO₂
FW: 431.3
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 273 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 1078 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 1078 in the solvent of choice, which should be purged with an inert gas. SR 1078 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SR 1078 in ethanol is approximately 15 mg/ml and approximately 10 mg/ml in DMSO and DMF.

SR 1078 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SR 1078 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SR 1078 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Retinoic acid receptor-related orphan receptors α and γ (ROR α and ROR γ) have been demonstrated to play important roles in the regulation of metabolism, circadian rhythms, immune function, and tumorigenesis. SR 1078 is a selective agonist of ROR α and ROR γ that stimulates ROR transcriptional activity in HEK293 cell reporter assays at concentrations as low as 2 μ M without effect at the related liver X receptors and farnesoid X receptors.¹ At 5 μ M, SR 1078 has been shown to stabilize p53 and to induce apoptosis in HepG2 cancer cells.²

References

1. Wang, Y., Kumar, N., Nuhant, P., *et al.* Identification of a synthetic agonist for the orphan nuclear receptors ROR α and ROR γ , SR1078. *ACS Chem. Biol.* **5**(11), 1029-1034 (2010).
2. Wang, Y., Solt, L.A., Kojetin, D.J., *et al.* Regulation of p53 stability and apoptosis by a ROR agonist. *PLoS One* **7**(4), 1-6 (2012).

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, 02/09/2024

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