

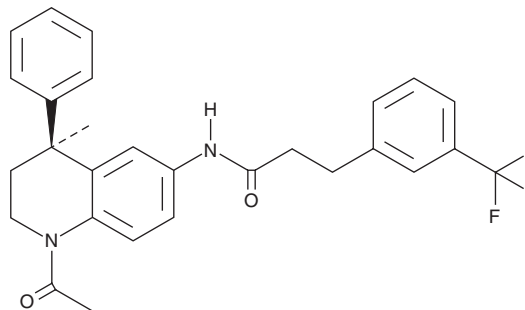
# PRODUCT INFORMATION



**ORG 274179-0**

Item No. 43557

**CAS Registry No.:** 1421683-12-6  
**Formal Name:** N-[(4S)-1-acetyl-1,2,3,4-tetrahydro-4-methyl-4-phenyl-6-quinolinyl]-3-(trifluoromethyl)benzenepropanamide  
**MF:** C<sub>28</sub>H<sub>27</sub>F<sub>3</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 480.5  
**Purity:** ≥98%  
**Supplied as:** A 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

ORG 274179-0 is supplied as a solid. A stock solution may be made by dissolving the ORG 274179-0 in the solvent of choice, which should be purged with an inert gas. ORG 274179-0 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

## Description

ORG 274179-0 is a thyroid-stimulating hormone receptor (TSHR) allosteric antagonist.<sup>1</sup> It inhibits TSHR activation induced by bovine TSH, human TSH, or the TSH-receptor-stimulating immunoglobulin M22 in reporter assays using CHO cells expressing the human receptor (IC<sub>50</sub>s = 30, 26, and 56 nM, respectively). ORG 274179-0 inhibits bovine TSH- or M22-induced phospholipase C (PLC) activation in CHO cells expressing human TSHR (IC<sub>50</sub>s = 5 and 4 nM, respectively).

## Reference

1. van Koppen, C.J., de Gooyer, M.E., Karstens, W.-J., *et al.* Mechanism of action of a nanomolar potent, allosteric antagonist of the thyroid-stimulating hormone receptor. *Br. J. Pharmacol.* **165(7)**, 2314-2324 (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, 08/12/2025

## 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