

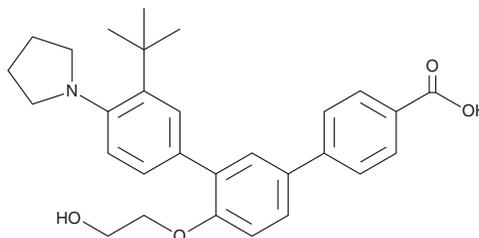
# PRODUCT INFORMATION



## Trifarotene

Item No. 45281

**CAS Registry No.:** 895542-09-3  
**Formal Name:** 3''-(1,1-dimethylethyl)-4'-(2-hydroxyethoxy)-4''-(1-pyrrolidinyl)-[1,1':3',1''-terphenyl]-4-carboxylic acid  
**Synonym:** CD5789  
**MF:** C<sub>29</sub>H<sub>33</sub>NO<sub>4</sub>  
**FW:** 459.6  
**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

Trifarotene is supplied as a solid. A stock solution may be made by dissolving the trifarotene in the solvent of choice, which should be purged with an inert gas. Trifarotene is soluble (≥10 mg/ml) in DMSO.

### Description

Trifarotene is a retinoic acid receptor  $\gamma$  (RAR $\gamma$ ) agonist.<sup>1</sup> It is selective for RAR $\gamma$  over RAR $\alpha$  and RAR $\beta$  in reporter assays using HeLa cells expressing chimeric RAR $\gamma$ , RAR $\alpha$ , or RAR $\beta$  receptors (EC<sub>50</sub>s = 7.7, 498, and 125 nM, respectively) and is inactive at the retinoid X receptor (RXR). Topical application of trifarotene (0.1% w/w) reduces ear edema induced by 12-O-tetradecanoylphorbol-1,3-acetate (TPA; Item No. 10008014) in a mouse model of epidermal inflammation. Formulations containing trifarotene have been used in the treatment of acne vulgaris.

### Reference

1. Aubert, J., Piwinica, D., Bertino, B., *et al.* Nonclinical and human pharmacology of the potent and selective topical retinoic acid receptor- $\gamma$  agonist trifarotene. *Br. J. Dermatol.* **179**(2), 442-456 (2018).

#### 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, 02/19/2026

#### 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