

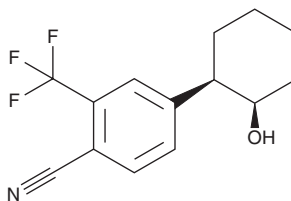
# PRODUCT INFORMATION



**PF-998425**

Item No. 31442

**CAS Registry No.:** 1076225-27-8  
**Formal Name:** 4-[(1R,2R)-2-hydroxycyclohexyl]-2-(trifluoromethyl)-benzonitrile  
**MF:** C<sub>14</sub>H<sub>14</sub>F<sub>3</sub>NO  
**FW:** 269.3  
**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

PF-998425 is supplied as a solid. A stock solution may be made by dissolving the PF-998425 in the solvent of choice, which should be purged with an inert gas. PF-998425 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

## Description

PF-998425 is a non-steroidal androgen receptor antagonist (IC<sub>50</sub> = 43 nM in a cell-based assay).<sup>1</sup> It binds to the androgen receptor (IC<sub>50</sub> = 37 nM in a radioligand binding assay) and is selective for the androgen receptor over the progesterone receptor (IC<sub>50</sub> = >10,000 nM). Topical administration of PF-998425 (3% twice per day) increases hair growth in a mouse model of hair growth, as well as reduces ear sebum in male Syrian golden hamsters (ED<sub>50</sub> = 0.4% twice per day).

## Reference

1. Li, J.J., Iula, D.M., Nguyen, M.N., *et al.* Rational design and synthesis of 4-[(1R,2R)-2-hydroxycyclohexyl]-2-(trifluoromethyl)benzonitrile (PF-998425), a novel, nonsteroidal androgen receptor antagonist devoid of phototoxicity for dermatological indications. *J. Med. Chem.* **51(21)**, 7010-7014 (2008).

### 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, 12/12/2022

## 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