

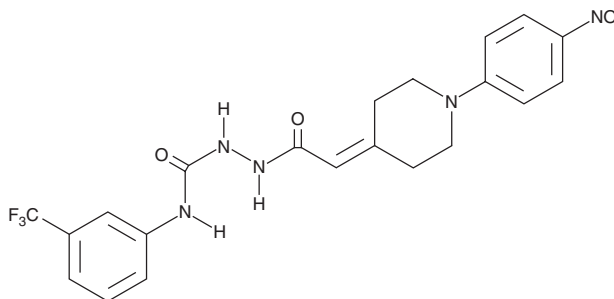
PRODUCT INFORMATION



Thyroid Hormone Receptor Antagonist (1-850)

Item No. 17248

CAS Registry No.: 251310-57-3
Formal Name: 2-[1-(4-nitrophenyl)-4-piperidinylidene]-acetic acid-2-[[[3-(trifluoromethyl)phenyl]amino]carbonyl]hydrazide
MF: C₂₁H₂₀F₃N₅O₄
FW: 463.4
Purity: ≥98%
UV/Vis.: λ_{max}: 253, 384 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

Thyroid hormone receptor antagonist (1-850) is supplied as a crystalline solid. A stock solution may be made by dissolving the thyroid hormone receptor antagonist (1-850) in the solvent of choice, which should be purged with an inert gas. Thyroid hormone receptor antagonist (1-850) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of thyroid hormone receptor antagonist (1-850) in ethanol is approximately 0.5 mg/ml and approximately 50 mg/ml in DMSO and DMF.

Thyroid hormone receptor antagonist (1-850) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, thyroid hormone receptor antagonist (1-850) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Thyroid hormone receptor antagonist (1-850) 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

Thyroid hormone receptor antagonist (1-850) competitively blocks T3 binding to both TR α and TR β (IC₅₀s = 1.5 μ M), demonstrating ~1,000-fold lower affinity for TR α in intact cells compared with T3.¹ This compound is specific, showing no effect on the activity of the related retinoic acid receptor α .¹

Reference

1. Schapira, M., Raaka, B.M., Das, S., *et al.* Discovery of diverse thyroid hormone receptor antagonists by high-throughput docking. *Proc. Natl. Acad. Sci. USA* **100**(12), 7354-7359 (2003).

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, 11/07/2022

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