

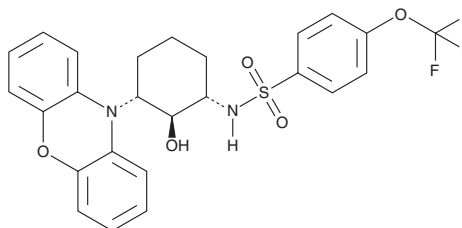
# PRODUCT INFORMATION



## (1S,2S,3R)-DT-061

Item No. 45615

**CAS Registry No.:** 1809427-20-0  
**Formal Name:** N-[(1S,2S,3R)-2-hydroxy-3-(10H-phenoxazin-10-yl)cyclohexyl]-4-(trifluoromethoxy)benzenesulfonamide  
**MF:** C<sub>25</sub>H<sub>23</sub>F<sub>3</sub>N<sub>2</sub>O<sub>5</sub>S  
**FW:** 520.5  
**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

(1S,2S,3R)-DT-061 is supplied as a solid. A stock solution may be made by dissolving the (1S,2S,3R)-DT-061 in the solvent of choice, which should be purged with an inert gas. (1S,2S,3R)-DT-061 is sparingly soluble (1-10 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

### Description

(1S,2S,3R)-DT-061 is the enantiomer of DT-061 (Item No. 42118).<sup>1</sup> It inhibits colony formation in NCI H1650 and A549 lung cancer cells (GI<sub>50</sub>s = 5 and 7.5 μM, respectively).

### Reference

1. Ohlmeyer, M. and Kastrinsky, D. Sulfonamides derived from tricyclic-2-aminocycloalkanol as anticancer agents. *Icahn School of Medicine at Mount Sinai* **US 9,937,186 B2** (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, 05/05/2026

#### 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