

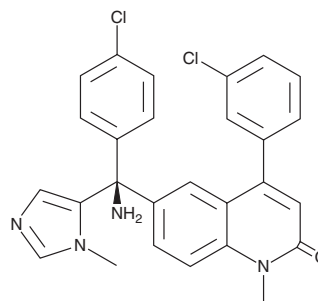
# PRODUCT INFORMATION



## (S)-Tipifarnib

Item No. 35515

**CAS Registry No.:** 192185-71-0  
**Formal Name:** 6-[(S)-amino(4-chlorophenyl)(1-methyl-1H-imidazol-5-yl)methyl]-4-(3-chlorophenyl)-1-methyl-2(1H)-quinolinone  
**Synonym:** (S)-(-)-R 115777  
**MF:** C<sub>27</sub>H<sub>22</sub>Cl<sub>2</sub>N<sub>4</sub>O  
**FW:** 489.4  
**Purity:** ≥98%  
**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

(S)-Tipifarnib is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-tipifarnib in the solvent of choice, which should be purged with an inert gas. (S)-Tipifarnib is soluble in organic solvents such as methanol and DMSO. The solubility of (S)-tipifarnib in methanol and DMSO is approximately 5 and 10 mg/ml, respectively.

### Description

(S)-Tipifarnib is an inactive enantiomer of tipifarnib (Item No. 11747).<sup>1</sup>

### Reference

1. Ren, P., Deng, X., and Mai, W. Synthesis of tipifarnib. *Kura Oncology, Inc.* **WO2019222565A1**, (2019).

#### 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, 10/25/2022

#### 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](http://WWW.CAYMANCHEM.COM)