

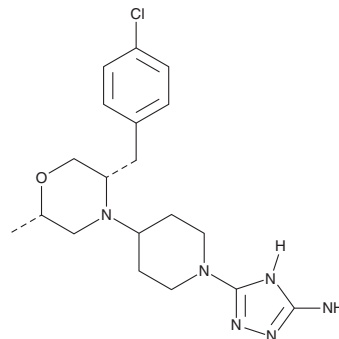
# PRODUCT INFORMATION



## OATD-01

Item No. 35193

**CAS Registry No.:** 2088453-21-6  
**Formal Name:** 5-[4-[(2S,5S)-5-[(4-chlorophenyl)methyl]-2-methyl-4-morpholinyl]-1-piperidinyl]-1H-1,2,4-triazol-3-amine  
**MF:** C<sub>19</sub>H<sub>27</sub>ClN<sub>6</sub>O  
**FW:** 390.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 225 nm  
**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

OATD-01 is supplied as a solid. A stock solution may be made by dissolving the OATD-01 in the solvent of choice, which should be purged with an inert gas. OATD-01 is soluble in ethanol.

### Description

OATD-01 is an inhibitor of chitotriosidase 1 (CHIT1) and acidic mammalian chitinase (CHIA), also known as AMCase (IC<sub>50</sub>s = 23 and 9 nM, respectively).<sup>1</sup> It is selective for CHIT1 and CHIA over a panel of 98 enzymes at 10 μM. OATD-01 (30 and 100 mg/kg) reduces the degree of lung fibrosis and plasma chitinolytic activity in a mouse model of pulmonary fibrosis induced by bleomycin (Item No. 13877).

### Reference

1. Koralweski, R., Dymek, B., Mazur, M., *et al.* Discovery of OATD-01, a first-in-class chitinase inhibitor as potential new therapeutics for idiopathic pulmonary fibrosis. *J. Med. Chem.* **63**(24), 15527-15540 (2020).

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