

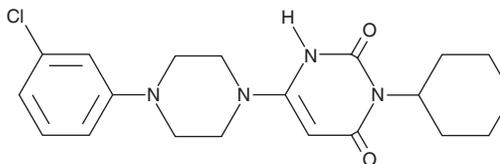
# PRODUCT INFORMATION



## ML-180

Item No. 15181

**CAS Registry No.:** 863588-32-3  
**Formal Name:** 6-[4-(3-chlorophenyl)-1-piperazinyl]-3-cyclohexyl-2,4(1H,3H)-pyrimidinedione  
**Synonyms:** CID-3238389, SR1848  
**MF:** C<sub>20</sub>H<sub>25</sub>ClN<sub>4</sub>O<sub>2</sub>  
**FW:** 388.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212, 274 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

ML-180 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-180 in the solvent of choice. ML-180 is soluble in DMSO at a concentration of approximately 10 mg/ml.

### Description

ML-180 is an inverse agonist of liver receptor homolog-1 (LRH-1, IC<sub>50</sub> = 3.7 μM) with maximum efficacy of 64% repression.<sup>1</sup> It is inactive at the related steroidogenic factor-1 transcriptional activator.<sup>1</sup> Through LRH-1, ML-180 alters the expression of haptoglobin and serum amyloid proteins A1 and A4, induces the death of estrogen-receptor negative MDA-MB-231 breast cancer cells, and inhibits the steroidogenic acute regulatory promoter (IC<sub>50</sub> = 2.05 μM).<sup>1</sup>

### Reference

1. Busby, S., Nuhant, P., Cameron, M., et al. Discovery of inverse agonists for the liver receptor homologue-1 (LRH1; NR5A2), in *Probe Reports from the NIH Molecular Libraries Program*,1 (2010).

#### 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, 11/28/2022

#### 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