

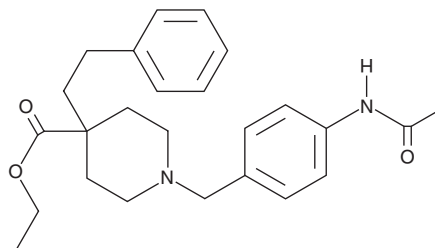
# PRODUCT INFORMATION



## ML-335

Item No. 15196

**CAS Registry No.:** 1069498-96-9  
**Formal Name:** 1-[[4-(acetylamino)phenyl]methyl]-4-(2-phenylethyl)-4-piperidinecarboxylic acid, ethyl ester  
**Synonyms:** CID-23723457, CYM 51010  
**MF:** C<sub>25</sub>H<sub>32</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 408.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 249 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-335 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-335 in the solvent of choice, which should be purged with an inert gas. ML-335 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

### Description

ML-335 is the first identified agonist for OPRM1-OPRD1 heterodimerization (EC<sub>50</sub> = 403 nM) that demonstrates selectivity over the individual OPRM1 and OPRD1 receptors and the serotonin HT5A receptor (EC<sub>50</sub>s = 14.9, 1.1, and >40 μM, respectively).<sup>1</sup>

### Reference

1. Pinello, C., Guerrero, M., Eberhart, C., et al. Characterization of an agonist probe for opioid receptor mu 1 (OPRM1)-opioid receptor delta 1 (OPRD1) heterodimerization. *Probe Reports from the NIH Molecular Libraries Program* 1-14 (2012).

#### 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, 12/16/2022

#### 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