

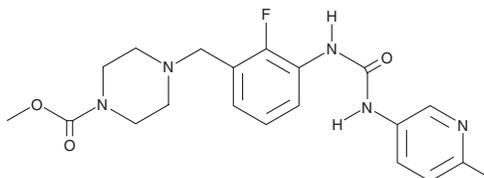
# PRODUCT INFORMATION



## Omecamtiv Mecarbil

Item No. 21074

**CAS Registry No.:** 873697-71-3  
**Formal Name:** 4-[[2-fluoro-3-[[[(6-methyl-3-pyridinyl)amino]carbonyl]amino]phenyl]methyl]-1-piperazinecarboxylic acid, methyl ester  
**Synonym:** CK-1827452  
**MF:** C<sub>20</sub>H<sub>24</sub>FN<sub>5</sub>O<sub>3</sub>  
**FW:** 401.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 256 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

Omecamtiv mecarbil is supplied as a crystalline solid. A stock solution may be made by dissolving the omecamtiv mecarbil in the solvent of choice. Omecamtiv mecarbil is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of omecamtiv mecarbil in these solvents is approximately 5, 10, and 30 mg/ml, respectively.

Omecamtiv mecarbil is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, omecamtiv mecarbil should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Omecamtiv mecarbil has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Omecamtiv mecarbil is a cardiac myosin ATPase activator with a 40% activation concentration (AC<sub>40</sub>) value of 0.58 μM for the bovine enzyme.<sup>1</sup> It increases fractional shortening in rats in a dose-dependent manner.

### Reference

1. Morgan, B.P., Muci, A., Lu, P.P., *et al.* Discovery of *omecamtiv mecarbil* the first, selective, small molecule activator of cardiac myosin. *ACS Med. Chem. Lett.* **1**(9), 472-477 (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, 12/22/2022

#### 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