

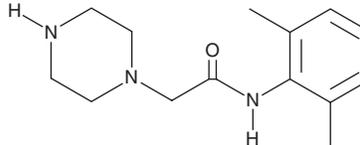
# PRODUCT INFORMATION



## CVT-2738

Item No. 35003

**CAS Registry No.:** 5294-61-1  
**Formal Name:** N-(2,6-dimethylphenyl)-1-piperazineacetamide  
**MF:** C<sub>14</sub>H<sub>21</sub>N<sub>3</sub>O  
**FW:** 247.3  
**Purity:** ≥98%  
**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

CVT-2738 is supplied as a solid. A stock solution may be made by dissolving the CVT-2738 in the solvent of choice, which should be purged with an inert gas. CVT-2738 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CVT-2738 in ethanol and DMSO is approximately 5 mg/ml and approximately 10 mg/ml in DMF.

CVT-2738 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CVT-2738 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CVT-2738 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

CVT-2738 is an active metabolite of the antianginal agent ranolazine (Item No. 15604).<sup>1</sup> It is formed from ranolazine by the cytochrome P450 (CYP) isoform CYP3A. CVT-2738 (100 mg/kg) inhibits increases in T-wave height and the RR interval in a mouse model of myocardial ischemia induced by isoprenaline (isoproterenol; Item No. 15592).<sup>2</sup>

### References

1. Wang, Y., Chen, X., Sun, Z., *et al.* Development and validation of a sensitive LC-MS/MS assay for simultaneous quantitation of ranolazine and its three metabolites in human plasma. *J. Chromatogr. B* **889-890**, 10-16 (2012).
2. Yao, Z., Gong, S., Guan, T., *et al.* Synthesis of ranolazine metabolites and their anti-myocardial ischemia activities. *Chem. Pharm. Bull. (Tokyo)* **57(11)**, 1218-1222 (2009).

#### 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, 12/09/2022

#### 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