

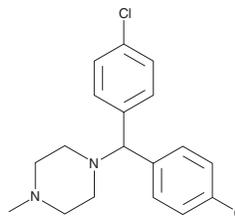
# PRODUCT INFORMATION



**CAY10704**

Item No. 19240

**CAS Registry No.:** 101784-44-5  
**Formal Name:** 1-[bis(4-chlorophenyl)methyl]-4-methyl-piperazine  
**MF:** C<sub>18</sub>H<sub>20</sub>Cl<sub>2</sub>N<sub>2</sub>  
**FW:** 335.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 233 nm  
**Supplied as:** A 5 mg/ml solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

CAY10704 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of CAY10704 in ethanol is 16 mg/ml and approximately 20 mg/ml in DMSO and DMF.

For maximum solubility in aqueous buffers, the DMF solution of CAY10704 should be diluted with the aqueous buffer of choice. CAY10704 has a solubility of 0.3 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

CAY10704 is a potent inhibitor of hepatitis C virus (HCV) infection (EC<sub>50</sub> = 17 nM) that displays low cytotoxicity of virally-infected human hepatoma Huh7.5.1 cells (CC<sub>50</sub> = 21.3 μM).<sup>1</sup> It displays good pharmacokinetics in mice when delivered intraperitoneally, with preferential liver distribution without significant hepatotoxicity.<sup>1</sup> CAY10704 is selective for HCV over dengue virus (EC<sub>50</sub> = 4.62 μM).<sup>1</sup>

## Reference

1. He, S., Xiao, J., Dulcey, A.E., *et al.* Discovery, optimization, and characterization of novel chlorcyclizine derivatives for the treatment of hepatitis C virus infection. *J. Med. Chem.* **59**(3), 841-853 (2016).

### 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, 01/05/2026

## 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