

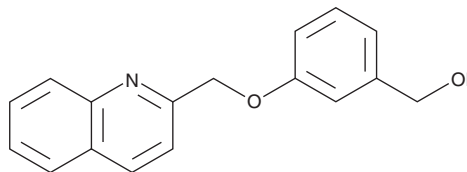
# PRODUCT INFORMATION



**CAY10789**

Item No. 35325

**CAS Registry No.:** 123226-28-8  
**Formal Name:** 3-(2-quinolinylmethoxy)-benzenemethanol  
**Synonym:** [3-(Quinolin-2-ylmethoxy)phenyl]methanol  
**MF:** C<sub>17</sub>H<sub>15</sub>NO<sub>2</sub>  
**FW:** 265.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 230 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

## Laboratory Procedures

CAY10789 is supplied as a solid. A stock solution may be made by dissolving the CAY10789 in the solvent of choice, which should be purged with an inert gas. CAY10789 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10789 in DMSO is approximately 5 mg/ml and approximately 11 mg/ml in ethanol and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CAY10789 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of CAY10789 in PBS (pH 7.2) is approximately 0.14 mg/ml. We do not recommend storing the aqueous solution for more than one day.

## Description

CAY10789 is an antagonist of the cysteinyl leukotriene 1 (CysLT<sub>1</sub>) receptor (IC<sub>50</sub> = 2.1 μM).<sup>1</sup>

## Reference

1. Fiorillo, B., Sepe, V., Conflitti, P., *et al.* Structural basis for developing multitarget compounds acting on cysteinyl leukotriene receptor 1 and G-protein-coupled bile acid receptor 1. *J. Med. Chem.* **64**(22), 16512-16529 (2021).

### 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, 06/01/2022

## 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