

# PRODUCT INFORMATION

## CM10

Item No. 36186

CAS Registry No.: 692269-09-3

Formal Name: 2-(2-propen-1-yl)-6-[[[1-(propyl-1H-benzimidazol-2-yl)amino]methyl]-phenol

MF:  $C_{20}H_{23}N_3O$

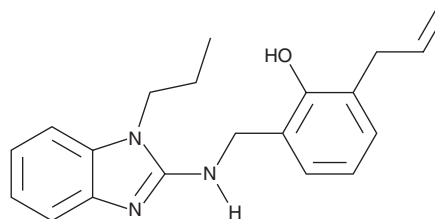
FW: 321.4

Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

CM10 is supplied as a solid. A stock solution may be made by dissolving the CM10 in the solvent of choice, which should be purged with an inert gas. CM10 is soluble in DMSO.

### Description

CM10 is a pan-aldehyde dehydrogenase 1 (ALDH1A) inhibitor ( $IC_{50}$ s = 1,700, 740, and 640 nM for human ALDH1A1, ALDH1A2, and ALDH1A3, respectively).<sup>1</sup> It is selective for ALDH1A isoforms over ALDH2, -3A1, -4A1, -5A1, -1B1, and -1L1 at 20  $\mu\text{M}$ . CM10 induces necroptosis in ovarian cancer stem cells.

### Reference

1. Chefetz, I., Grimley, E., Yang, K., *et al.* A Pan-ALDH1A inhibitor induces necroptosis in ovarian cancer stem-like cells. *Cell Rep.* **26(11)**, 3061-3075 (2019).

#### 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, 12/06/2022

#### 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