

# PRODUCT INFORMATION

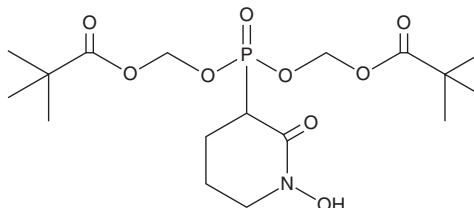


## POMHEX

Item No. 33390

**CAS Registry No.:** 2004714-34-3  
**Formal Name:** 2,2-dimethyl-propanoic acid, 1,1'-[[[(1-hydroxy-2-oxo-3-piperidiny)phosphinylidene]bis(oxymethylene)] ester

**MF:** C<sub>17</sub>H<sub>30</sub>NO<sub>9</sub>P  
**FW:** 423.4  
**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

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

### Description

POMHEX is a prodrug form of the  $\beta$ -enolase, also known as enolase-3, inhibitor HEX.<sup>1</sup> It is converted to HEX by intracellular carboxylesterases and selectively reduces the cell density of *ENO1*<sup>-/-</sup> D-423MG glioblastoma cells (IC<sub>50</sub> = 38 nM) over *ENO1*-overexpressing D-423MG cells and LN-319 astrocytoma cells expressing wild-type *ENO1* (IC<sub>50</sub>s = 604 and 2,148 nM, respectively). POMHEX decreases cell density in normoxic and hypoxic growth conditions in *ENO1*<sup>-/-</sup> D-423MG cells (IC<sub>50</sub>s = 79 and 201 nM, respectively).

### Reference

1. Yan, V.C., Yang, K.L., Ballato, E.S., *et al.* Bioreducible phosphonoamidate pro-drug inhibitor of enolase: Proof of concept study. *ACS Med. Chem. Lett.* **11**(7), 1484-1489 (2020).

#### 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, 08/04/2023

#### 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