

# PRODUCT INFORMATION

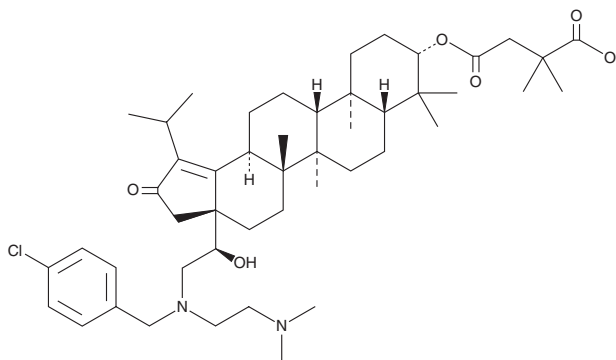


**GSK2838232**

Item No. 21390

**CAS Registry No.:** 1443460-91-0  
**Formal Name:** (3 $\beta$ )-3-(3-carboxy-3-methyl-1-oxobutoxy)-17-[(1R)-2-[[[4-chlorophenyl)methyl][2-(dimethylamino)ethyl]amino]-1-hydroxyethyl]-28-norlup-18-en-21-one

**MF:** C<sub>48</sub>H<sub>73</sub>ClN<sub>2</sub>O<sub>6</sub>  
**FW:** 809.6  
**Purity:**  $\geq 95\%$   
**UV/Vis.:**  $\lambda_{\text{max}}$ : 222, 247 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



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

## Laboratory Procedures

GSK2838232 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2838232 in the solvent of choice, which should be purged with an inert gas. GSK2838232 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of GSK2838232 in these solvents is approximately 1 mg/ml.

GSK2838232 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK2838232 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. GSK2838232 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

GSK2838232 is a second generation inhibitor of HIV-1 maturation ( $IC_{90} = 6.4$  nM).<sup>1</sup> It is active against a variety of HIV isolates resistant to bevirimat (PA-457; Item No. 21339).<sup>2</sup>

## References

1. Johnson, M., Jewell, R.C., Peppercorn, A., *et al.* The safety, tolerability, and pharmacokinetic profile of GSK2838232, a novel 2nd generation HIV maturation inhibitor, as assessed in healthy subjects. *Pharmacol. Res. Perspect.* **6**(4), e00408 (2018).
2. Cihlar, T. and Fordyce, M. Current status and prospects of HIV treatment. *Curr. Opin. Virol.* **18**, 50-56 (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, 11/30/2022

## 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](http://WWW.CAYMANCHEM.COM)