

# PRODUCT INFORMATION

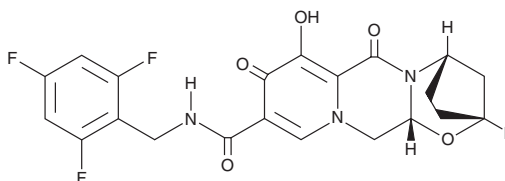


## Bictegravir

Item No. 26532

**CAS Registry No.:** 1611493-60-7  
**Formal Name:** (2R,5S,13aR)-2,3,4,5,7,9,13,13a-octahydro-8-hydroxy-7,9-dioxo-N-[(2,4,6-trifluorophenyl)methyl]-2,5-methanopyrido[1',2':4,5]pyrazino[2,1-b][1,3]oxazepine-10-carboxamide

**Synonym:** GS-9883  
**MF:** C<sub>21</sub>H<sub>18</sub>F<sub>3</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 449.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 248, 260, 322 nm  
**Supplied as:** A crystalline 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

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

### Description

Bictegravir is an inhibitor of HIV-1 integrase (IC<sub>50</sub> = 7.5 nM for strand transfer activity).<sup>1</sup> It has antiviral activity against clinical isolates of HIV-1 in human peripheral blood mononuclear cells (PBMCs; EC<sub>50</sub>s = 0.04-1.7 nM). Bictegravir also inhibits HIV-1 viral infection in MT-2 and MT-4 cells, CD4<sup>+</sup> T cells, and macrophages (EC<sub>50</sub>s = 1.5, 2.5, 1.5, and 6.6 nM, respectively) without exhibiting cytotoxicity (CC<sub>50</sub>s = 10.3, 3.7, 13, and 29.8 μM, respectively). Formulations containing bictegravir in combination with emtricitabine and tenofovir alafenamide have been used in the treatment of HIV-1 infections.

### Reference

1. Tsiang, M., Jones, G.S., Goldsmith, J., *et al.* Antiviral activity of bictegravir (GS-9883), a novel potent HIV-1 integrase strand transfer inhibitor with an improved resistance profile. *Antimicrob. Agents Chemother.* **60**(12), 7086-7097 (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, 10/18/2023

#### 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