

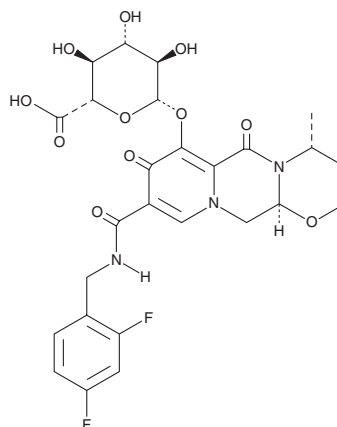
PRODUCT INFORMATION



Dolutegravir O-β-D-Glucuronide

Item No. 34509

CAS Registry No.: 1485692-21-4
Formal Name: (4R,12aS)-9-[[[(2,4-difluorophenyl)methyl]amino]carbonyl]-3,4,6,8,12,12a-hexahydro-4-methyl-6,8-dioxo-2H-pyrido[1',2':4,5]pyrazino[2,1-b][1,3]oxazin-7-yl β-D-glucopyranosiduronic acid
Synonyms: Dolutegravir Glucuronide, DTG Glucuronide
MF: C₂₆H₂₇F₂N₃O₁₁
FW: 595.5
Purity: ≥95%
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

Dolutegravir O-β-D-glucuronide is supplied as a solid. Dolutegravir O-β-D-glucuronide is slightly soluble in water (sonicated). We do not recommend storing the aqueous solution for more than one day.

Description

Dolutegravir O-β-D-glucuronide is a metabolite of the HIV integrase inhibitor dolutegravir (Item No. 22191).¹ It is formed from dolutegravir primarily by the UDP-glucuronosyltransferase (UGT) isoform UGT1A1 *in vivo* but is also metabolized by UGT1A9 in human liver and kidney microsomes and UGT1A3 in human intestinal microsomes.^{1,2}

References

1. Liu, S.N., Lu, J.B., Watson, C.J.W., *et al.* Mechanistic assessment of extrahepatic contributions to glucuronidation of integrase strand transfer inhibitors. *Drug Metab. Dispos.* **47(5)**, 535-544 (2019).
2. Reese, M.J., Savina, P.M., Generaux, G.T., *et al.* In vitro investigations into the roles of drug transporters and metabolizing enzymes in the disposition and drug interactions of dolutegravir, a HIV integrase inhibitor. *Drug Metab. Dispos.* **41(2)**, 353-361 (2013).

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, 10/21/2022

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