

PRODUCT INFORMATION

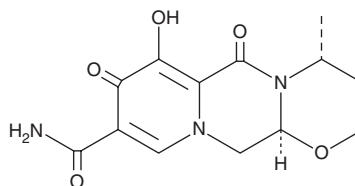


Dolutegravir M1

Item No. 34508

CAS Registry No.: 1485081-26-2
Formal Name: 3,4R,6,8,12,12aS-hexahydro-7-hydroxy-4-methyl-6,8-dioxo-2H-pyrido[1',2':4,5]pyrazino[2,1-b][1,3]oxazine-9-carboxamide

MF: C₁₃H₁₅N₃O₅
FW: 293.3
Purity: ≥95%
UV/Vis.: λ_{max}: 257, 323 nm
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 M1 is supplied as a solid. A stock solution may be made by dissolving the dolutegravir M1 in the solvent of choice, which should be purged with an inert gas. Dolutegravir M1 is slightly soluble in methanol.

Description

Dolutegravir M1 is a metabolite of the HIV integrase inhibitor dolutegravir (Item No. 22191).¹ It is formed primarily by the cytochrome P450 (CYP) isoforms CYP1A1 and CYP1B1.

Reference

1. Zhu, J., Wang, P., Li, F., *et al.* CYP1A1 and 1B1-mediated metabolic pathways of dolutegravir, an HIV integrase inhibitor. *Biochem. Pharmacol.* **158**, 174-184 (2018).

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, 10/28/2022

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