

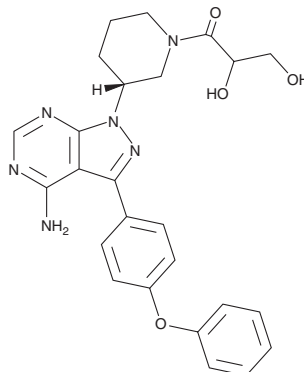
PRODUCT INFORMATION



PCI 45227

Item No. 33560

CAS Registry No.: 1654820-87-7
Formal Name: 1-[(3R)-3-[4-amino-3-(4-phenoxyphenyl)-1H-pyrazolo[3,4-d]pyrimidin-1-yl]-1-piperidinyl]-2,3-dihydroxy-1-propanone
Synonym: Dihydrodiol Ibrutinib
MF: C₂₅H₂₆N₆O₄
FW: 474.5
Purity: ≥95% (mixture of diastereomers)
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

PCI 45227 is supplied as a solid. A stock solution may be made by dissolving the PCI 45227 in the solvent of choice, which should be purged with an inert gas. PCI 45227 is soluble in methanol.

Description

PCI 45227 is an active metabolite of the Bruton's tyrosine kinase inhibitor ibrutinib (Item No. 16274).¹ PCI 45227 is formed from ibrutinib by the cytochrome P450 (CYP) isoform CYP3A.

Reference

1. Veeraraghavan, S., Viswanadha, S., Thappali, S., *et al.* Simultaneous quantification of lenalidomide, ibrutinib and its active metabolite PCI-45227 in rat plasma by LC-MS/MS: Application to a pharmacokinetic study. *J. Pharm. Biomed. Anal.* **107**, 151-158 (2015).

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, 11/09/2022

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