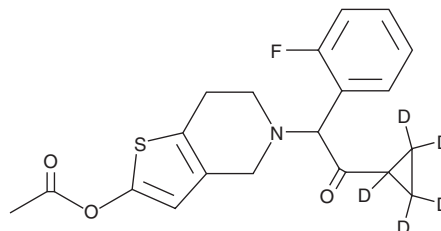


PRODUCT INFORMATION



Prasugrel-d₅ Item No. 28160

CAS Registry No.: 1127252-92-9
Formal Name: 2-[2-(acetyloxy)-6,7-dihydrothieno[3,2-c]pyridin-5(4H)-yl]-1-(cyclopropyl-1,2,2,3,3-d₅)-2-(2-fluorophenyl)-ethanone
MF: C₂₀H₁₅D₅FNO₃S
FW: 378.5
Chemical Purity: ≥95% (Prasugrel)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
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

Prasugrel-d₅ is intended for use as an internal standard for the quantification of prasugrel (Item No. 14278) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Prasugrel-d₅ is supplied as a solid. A stock solution may be made by dissolving the prasugrel-d₅ in the solvent of choice, which should be purged with an inert gas. Prasugrel-d₅ is slightly soluble in methanol and chloroform.

Description

Prasugrel is a prodrug form of the platelet purinergic P2Y₁₂ receptor antagonist R-99224.¹ Prasugrel (0.3 and 3 mg/kg) inhibits *ex vivo* washed platelet aggregation in rat platelet rich-plasma.² *In vivo*, prasugrel prevents thrombus formation (ED₅₀ = 0.68 mg/kg) and increases tail bleeding time in rats. Formulations containing prasugrel have been used in the prevention of blood clots.

References

1. Sugidachi, A., Asai, F., Yoneda, K., *et al.* Antiplatelet action of R-99224, an active metabolite of a novel thienopyridine-type G_i-linked P2T antagonist, CS-747. *Br. J. Pharmacol.* **132(1)**, 47-54 (2001).
2. Sugidachi, A., Asai, F., Ogawa, T., *et al.* The *in vivo* pharmacological profile of CS-747, a novel antiplatelet agent with platelet ADP receptor antagonist properties. *Br. J. Pharmacol.* **129(7)**, 1439-1446 (2000).

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, 12/19/2022

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