

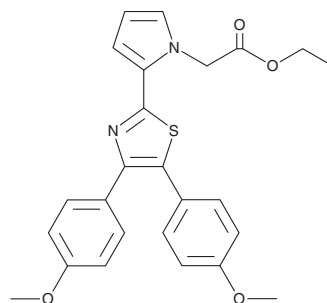
PRODUCT INFORMATION



Pamicogrel

Item No. 33514

CAS Registry No.: 101001-34-7
Formal Name: 2-[4,5-bis(4-methoxyphenyl)-2-thiazolyl]-1H-pyrrole-1-acetic acid, ethyl ester
Synonym: KBT-3022
MF: C₂₅H₂₄N₂O₄S
FW: 448.5
Purity: ≥98%
UV/Vis.: λ_{max}: 289, 347 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

Pamicogrel is supplied as a crystalline solid. A stock solution may be made by dissolving the pamicogrel in the solvent of choice, which should be purged with an inert gas. Pamicogrel is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of pamicogrel in these solvents is approximately 5 and 10 mg/ml, respectively.

Pamicogrel is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pamicogrel should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Pamicogrel has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Pamicogrel is a COX inhibitor and an antiplatelet agent.¹⁻³ It inhibits platelet aggregation induced by arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) or collagen in washed isolated rabbit platelets (IC₅₀s = 1.5 and 0.43 μM, respectively).² Pamicogrel (1 mg/kg) reduces thrombus weight in a guinea pig model of arteriovenous shunt thrombosis.³

References

1. Light, P. Pamicogrel (Kanebo). *IDrugs* **2(9)**, 938-942 (1999).
2. Yokota, K., Yamamoto, N., Morimoto, Y., *et al.* Effect of KBT-3022, a new diphenylthiazole derivative, on platelet functions. *J. Pharm. Pharmacol.* **47(9)**, 768-774 (1995).
3. Yokota, K., Yamashita, A., and Oda, M. Anti-thrombotic activity of KBT-3022 in experimental models of thrombosis. *Jpn. J. Pharmacol.* **68(2)**, 201-206 (1995).

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, 11/23/2022

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