

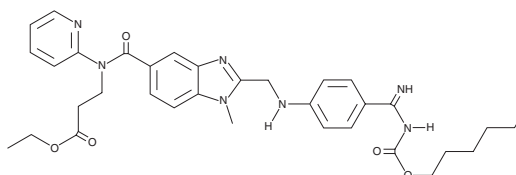
PRODUCT INFORMATION



Dabigatran etexilate

Item No. 17131

CAS Registry No.: 211915-06-9
Formal Name: N-[[2-[[[4-[[[(hexloxy)carbonyl]amino]iminomethyl]phenyl]amino]methyl]-1-methyl-1H-benzimidazol-5-yl]carbonyl]-N-2-pyridinyl-b-alanine, ethyl ester
Synonym: BIBR 1048
MF: C₃₄H₄₁N₇O₅
FW: 627.7
Purity: ≥95%
UV/Vis.: λ_{max}: 209, 227, 318 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

Dabigatran etexilate is supplied as a crystalline solid. A stock solution may be made by dissolving the dabigatran etexilate in the solvent of choice, which should be purged with an inert gas. Dabigatran etexilate is soluble in the organic solvent DMSO at a concentration of approximately 0.5 mg/ml at 25°C.

Description

Dabigatran etexilate is a prodrug form of the thrombin inhibitor dabigatran (Item No. 17133).¹ It inhibits thrombin-induced platelet aggregation in isolated rabbit platelets (IC₅₀ = 0.337 μM) and inhibits thrombus formation in a rat model of venous thrombosis (ED₅₀ = 7.8 μmol/kg). Formulations containing dabigatran etexilate have been used to reduce the risk of stroke and systemic embolism in patients with non-valvular atrial fibrillation and in the treatment and prevention of deep vein thrombosis (DVT) and pulmonary embolism.

Reference

1. Yang, X.-Z., Diao, X.-J., Yang, W.-H., *et al.* Design, synthesis and antithrombotic evaluation of novel dabigatran prodrugs containing methyl ferulate. *Bioorg. Med. Chem. Lett.* **23(7)**, 2089-2092 (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, 11/11/2022

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