

PRODUCT INFORMATION



Dabigatran Acyl-β-D-Glucuronide (trifluoroacetate salt)

Item No. 39046

Formal Name: N-[[2-[[[4-(aminoiminomethyl)phenyl]amino)methyl]-1-methyl-1H-benzimidazol-5-yl]carbonyl]-N-2-pyridinyl-β-alanine, β-D-glucopyranuronosyl ester, trifluoroacetate salt

Synonym: Dabigatran β-1-O-Acyl Glucuronide

MF: C₃₁H₃₃N₇O₉ • XCF₃COOH

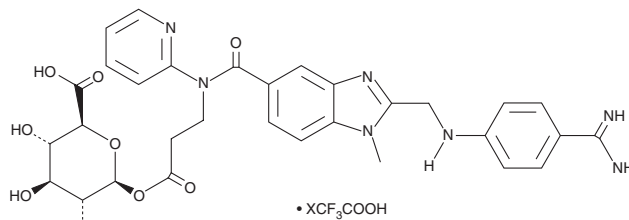
FW: 647.6

Purity: ≥75%

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

Dabigatran acyl-β-D-glucuronide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the dabigatran acyl-β-D-glucuronide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Dabigatran acyl-β-D-glucuronide (trifluoroacetate salt) is slightly soluble in methanol and DMSO.

Description

Dabigatran acyl-β-D-glucuronide is an active metabolite of the thrombin inhibitor dabigatran (Item No. 17133).^{1,2} It is formed from dabigatran primarily by the UDP-glucuronosyltransferase (UGT) isoform UGT2B15. Dabigatran acyl-β-D-glucuronide (0.46 μM) increases activated partial thromboplastin time (aPTT) in isolated human platelet-poor plasma.²

References

1. Eisert, W.G., Huel, N., Stangier, J., *et al.* Dabigatran: An oral novel potent reversible nonpeptide inhibitor of thrombin. *Arterioscler. Thromb. Vasc. Biol.* **30(10)**, 1885-1889 (2010).
2. Ebner, T., Wagner, K., and Wiene, W. Dabigatran acylglucuronide, the major human metabolite of dabigatran: In vitro formation, stability, and pharmacological activity. *Drug. Metab. Dispos.* **38(9)**, 1567-1575 (2010).

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, 06/27/2023

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