

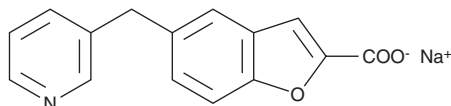
PRODUCT INFORMATION



Furegrelate (sodium salt)

Item No. 70540

CAS Registry No.: 85666-17-7
Formal Name: 5-(3-pyridinylmethyl)-2-benzofurancarboxylic acid, sodium salt
Synonym: U-63557A
MF: C₁₅H₁₀NO₃ • Na
FW: 275.2
Purity: ≥99%
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

Furegrelate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the furegrelate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Furegrelate (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of furegrelate (sodium salt) in these solvents is approximately 14.5, 21, and 25 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of furegrelate (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of furegrelate (sodium salt) in PBS (pH 7.2) is approximately 16.5. We do not recommend storing the aqueous solution for more than one day.

Description

Furegrelate is a potent inhibitor of thromboxane synthase with little effect on other enzymes essential for arachidonate metabolism. The IC₅₀ value is 15 nM for human platelet microsomal thromboxane synthase.^{1,2}

References

1. Johnson, R.J., Nidy, E.G., Aiken, J.W., *et al.* Thromboxane A₂ synthase inhibitors. 5-(3-pyridylmethyl) benzofuran-2-carboxylic acids. *J. Med. Chem.* **29**, 1461-1468 (1986).
2. Gorman, R.R., Johnson, R.A., Spilman, C.H., *et al.* Inhibition of platelet thromboxane A₂ synthase activity by sodium 5-(3'-pyridinylmethyl)benzofuran-2-carboxylate. *Prostaglandins* **26**, 325-342 (1983).

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, 04/05/2024

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