

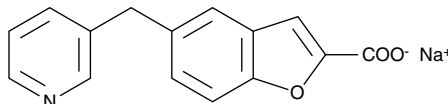
Product Information



Furegrelate (sodium salt)

Item No. 70540

CAS Registry No: 85666-17-7
Formal Name: 5-(3-pyridinylmethyl)-2-benzofuran-2-carboxylic acid, sodium salt
Synonym: U-63557A
MF: C₁₅H₁₀NO₃ • Na
FW: 275.2
Purity: ≥99%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
Melting Point: 347-349°C



Laboratory Procedures

For long term storage, we suggest that furegrelate (sodium salt) be stored as supplied at -20°C. It should be stable for at least one year.

Before use, a stock solution should be prepared by dissolving furegrelate (sodium salt) in water. The solubility of furegrelate (sodium salt) in water is approximately 15 mg/ml. Stock solutions will be stable for at least a month if stored at -20°C in an oxygen-free environment.

Furegrelate is a potent inhibitor of thromboxane (TX) synthase, and has no effect on other enzymes in arachidonate metabolism.¹ At a dose of 3.0 mg/kg in rhesus monkeys, furegrelate inhibited TX synthase by 80% for at least 12 hours.¹ *In vitro* experiments give an ED₅₀ range of 2-800 ng/ml, depending on the source of TX synthase.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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/70540

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 08/08/2012

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com