

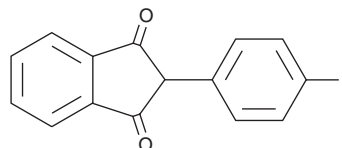
PRODUCT INFORMATION



Fluindione

Item No. 34966

CAS Registry No.: 957-56-2
Formal Name: 2-(4-fluorophenyl)-1H-indene-1,3(2H)-dione
MF: $C_{15}H_9FO_2$
FW: 240.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 223, 267 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Fluindione is supplied as a solid. A stock solution may be made by dissolving the fluindione in the solvent of choice, which should be purged with an inert gas. Fluindione is soluble in ethanol and DMSO. The solubility of fluindione in these solvents is approximately 1 mg/ml.

Description

Fluindione is a vitamin K antagonist.² It inhibits vitamin K epoxide reductase (VKOR; $IC_{50} = 6.6$ nM), including VKOR containing certain substitution mutations. Fluindione inhibits platelet and fibrin deposition, as well as prevents β -thromboglobulin release and thrombin-antithrombin complex formation, in an ex vivo model of tissue factor-induced thrombus formation using isolated human whole blood.¹ Fluindione has previously been used in the treatment of emboligenic heart disease, deep vein thrombosis, pulmonary embolism, and in the prevention of thromboembolic complications in patients with heart disease.

References

1. Bossavy, J.P., Sakariassen, K.S., Thalamas, C., *et al.* Antithrombotic efficacy of the vitamin K antagonist fluindione in a human ex vivo model of arterial thrombosis: Effect of anticoagulation level and combination therapy with aspirin. *Arterioscler. Thromb. Vasc. Biol.* **19**(9), 2269-2275 (1999).
2. Chen, X., Jin, D.-Y., Stafford, D.W., *et al.* Evaluation of oral anticoagulants with vitamin K epoxide reductase in its native milieu. *Blood* **132**(18), 1974-1984 (2018).

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, 11/14/2022

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