

PRODUCT INFORMATION



Clanfener

Item No. 43105

CAS Registry No.: 51213-99-1

Formal Name: N-[[[4-chlorophenyl]amino]carbonyl]-2-(dimethylamino)-6-fluoro-benzamide

Synonym: NSC 603108

MF: $C_{16}H_{15}ClFN_3O_2$

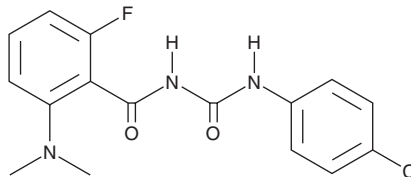
FW: 335.8

Purity: $\geq 98\%$

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

Clanfener is supplied as a solid. A stock solution may be made by dissolving the clanfener in the solvent of choice, which should be purged with an inert gas. Clanfener is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Clanfener is an anticancer agent.¹ It induces cell death in B16/F10 melanoma cells with a half-maximal inhibitory dose (ID_{50}) value of 1 $\mu\text{g/ml}$. Clanfener (800 mg/kg) also increases the number of bone marrow cells, mature and immature neutrophils, and decreases the number of proerythroblasts and normoblast erythrocytes in mice after 42 hours.²

References

1. Hofs, H.P. and McVie, J.G. Effects of benzoylphenyl ureas on growth of B16 melanoma cells *in vitro* and *in vivo*. *Invest. New Drugs* **9**(3), 227-232 (1991).
2. Jenkins, V.K., Juneja, H.S., Ives, K., *et al.* Effects of diflubenzuron and clanfener on mouse bone marrow cells. *Invest. New Drugs* **11**(4), 279-289 (1993).

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, 05/08/2025

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