

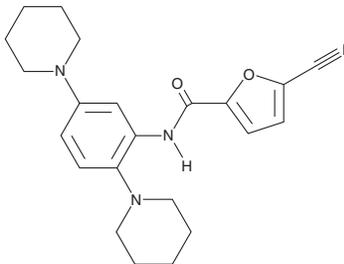
PRODUCT INFORMATION



cFMS Receptor Tyrosine Kinase Inhibitor

Item No. 44282

CAS Registry No.: 959626-45-0
Formal Name: 5-cyano-N-(2,5-di-1-piperidinylphenyl)-2-furancarboxamide
MF: C₂₂H₂₆N₄O₂
FW: 378.5
Purity: ≥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

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

Description

cFMS receptor tyrosine kinase inhibitor is an inhibitor of CSF-1 receptor tyrosine kinase (FMS; IC₅₀ = 0.017 μM).¹

Reference

1. Patch, R.J., Brandt, B.M., Asgari, D., *et al.* Potent 2'-aminoanilide inhibitors of cFMS as potential anti-inflammatory agents. *Bioorg. Med. Chem. Lett.* **17(22)**, 6070-6074 (2007).

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, 09/10/2025

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