

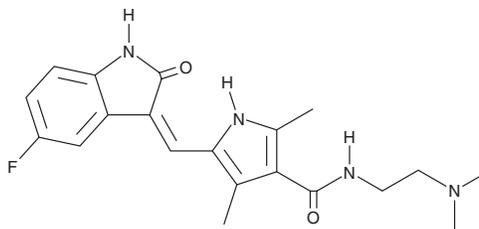
PRODUCT INFORMATION



N,N-dimethyl Sunitinib

Item No. 45148

CAS Registry No.: 326914-17-4
Formal Name: N-[2-(dimethylamino)ethyl]-5-[(Z)-(5-fluoro-1,2-dihydro-2-oxo-3H-indol-3-ylidene)methyl]-2,4-dimethyl-1H-pyrrole-3-carboxamide
Synonym: SU 11657
MF: C₂₀H₂₃FN₄O₂
FW: 370.4
Purity: ≥95%
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

N,N-dimethyl Sunitinib is supplied as a solid. A stock solution may be made by dissolving the N,N-dimethyl sunitinib in the solvent of choice, which should be purged with an inert gas. N,N-dimethyl Sunitinib is sparingly soluble (1-10 mg/ml) in DMSO.

Description

N,N-dimethyl Sunitinib is a multi-kinase inhibitor and derivative of sunitinib (Item No. 13159).¹ It inhibits VEGFR1, VEGFR2, PDGFR, and c-Kit (IC₅₀s = 0.013, 0.017, 0.005, and 0.04 μM, respectively). Oral administration of N,N-dimethyl sunitinib (40 mg/kg) reduces the degradation of cartilage proteoglycan and the severity of arthritis in the K/BxN mouse model of rheumatoid arthritis. N,N-dimethyl Sunitinib also reduces intratumoral angiogenesis and tumor growth in SK-N-AS, MYCN-amplified IMR-32, and SH-SY5Y neuroblastoma mouse xenograft models.²

References

1. De Bandt, M., Ben Mahdi, M.H., Ollivier, V., *et al.* Blockade of vascular endothelial growth factor receptor I (VEGF-RI), but not VEGF-RII, suppresses joint destruction in the K/BxN model of rheumatoid arthritis. *J. Immunol.* **171(9)**, 4853-4859 (2003).
2. Bäckman, U. and Christofferson, R. The selective class III/V receptor tyrosine kinase inhibitor SU11657 inhibits tumor growth and angiogenesis in experimental neuroblastomas grown in mice. *Pediatr. Res.* **57(5 Pt 1)**, 690-695 (2005).

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, 03/03/2026

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