

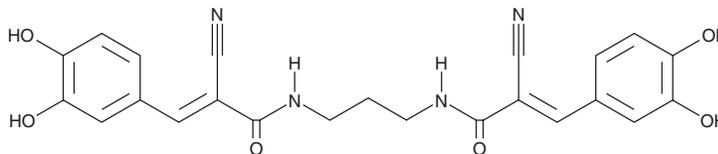
PRODUCT INFORMATION



AG-537

Item No. 45255

CAS Registry No.: 140674-77-7
Formal Name: (2E,2'E)-N,N'-1,3-propanediylbis[2-cyano-3-(3,4-dihydroxyphenyl)-2-propenamamide]
Synonyms: Tyrphostin AG 537, Tyrphostin 537
MF: C₂₃H₂₀N₄O₆
FW: 448.4
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

AG-537 is supplied as a solid. A stock solution may be made by dissolving the AG-537 in the solvent of choice, which should be purged with an inert gas. AG-537 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

Description

AG-537 is a dimeric tyrphostin inhibitor of EGFR (IC₅₀ = 0.4 μM).¹ It is selective for EGFR over the Bcr-Abl fusion proteins p210 and p185 and wild-type c-Abl (IC₅₀s = 44, >100, and >100 μM, respectively, for the human proteins). AG-537 inhibits EGFR autophosphorylation (IC₅₀ = 3.3 μM) and EGF-dependent proliferation of HER14 fibroblasts (IC₅₀ = 3 μM).² It is also an inhibitor of *T. brucei* RNA-editing ligase 1 (REL1; IC₅₀ = 0.25 μM) and *M. tuberculosis* 4'-phosphopantetheinyl transferase (PptT; EC₅₀ = 5.9 μM in an apo-blue pigment synthetase A (BpsA) activation assay).^{3,4}

References

1. Anafi, M., Gazit, A., Gilon, C., *et al.* Selective interactions of transforming and normal abl proteins with ATP, tyrosine-copolymer substrates, and tyrphostins. *J. Biol. Chem.* **267**(8), 4518-4523 (1992).
2. Gazit, A., Oshero, N., Gilon, C., *et al.* Tyrphostins. 6. Dimeric benzylidenemalononitrile tyrphostins: Potent inhibitors of EGF receptor tyrosine kinase in vitro. *J. Med. Chem.* **39**, 4905-4911 (1996).
3. Zimmermann, S., Hall, L., Riley, S., *et al.* A novel high-throughput activity assay for the *Trypanosoma brucei* editosome enzyme REL1 and other RNA ligases. *Nucleic Acids Res.* **44**(3), e24 (2016).
4. Brown, A.S., Owen, J.G., Jung, J., *et al.* Inhibition of indigoidine synthesis as a high-throughput colourimetric screen for antibiotics targeting the essential *Mycobacterium tuberculosis* phosphopantetheinyl transferase PptT. *Pharmaceuticals* **13**(7), 1066 (2021).

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, 02/24/2026

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