

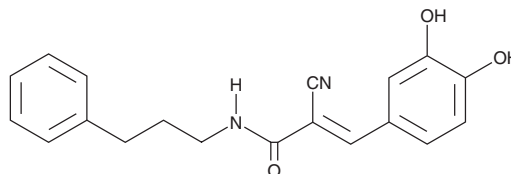
PRODUCT INFORMATION



AG-555

Item No. 23247

CAS Registry No.: 133550-34-2
Formal Name: (2E)-2-cyano-3-(3,4-dihydroxyphenyl)-N-(3-phenylpropyl)-2-propenamide
Synonyms: Tyrphostin AG-555, Tyrphostin B46
MF: C₁₉H₁₈N₂O₃
FW: 322.4
Purity: ≥95%
UV/Vis.: λ_{max}: 257, 362 nm
Supplied as: A crystalline 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-555 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-555 in the solvent of choice, which should be purged with an inert gas. AG-555 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AG-555 in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

AG-555 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AG-555 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AG-555 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AG-555 is a tyrphostin inhibitor of the EGF receptor (EGFR; IC₅₀ = 0.7 μM).¹ It selectively inhibits EGFR over ErbB2 (IC₅₀ = 35 μM). AG-555 inhibits EGF-dependent growth of HER 14 cells (IC₅₀ = 2.5 μM) as well as the growth of psoriatic keratinocytes isolated from patients with psoriasis when used at concentrations ranging from 1 to 50 μM.^{1,2} It also inhibits Moloney murine leukemia virus (Mo-MuLV) reverse transcriptase activity (IC₅₀ = 10.8 μM) without affecting Mo-MuLV-infected NIH3T3 cell growth (IC₅₀ = 210 μM).³

References

1. Gazit, A., Oshero, N., Posner, I., *et al.* Tyrphostins. 2. Heterocyclic and α-substituted benzylidenemalononitrile tyrphostins as potent inhibitors of EGF receptor and ErbB2/neu tyrosine kinases. *J. Med. Chem.* **34**(6), 1896-1907 (1991).
2. Ben-Bassat, H., Vardi, D.V., Gazit, A., *et al.* Tyrphostins suppress the growth of psoriatic keratinocytes. *Exp. Dermatol.* **4**(2), 82-88 (1995).
3. Esther, A., Iftach, S., and Esther, P. Inhibition of Moloney murine leukemia virus replication by tyrphostins, tyrosine kinase inhibitors. *FEBS Lett.* **341**(1), 99-103 (1994).

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, 11/17/2022

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