

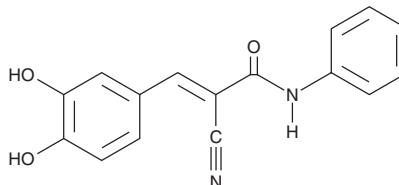
PRODUCT INFORMATION



AG-494

Item No. 10010242

CAS Registry No.: 133550-35-3
Formal Name: 2E-cyano-3-(3,4-dihydroxyphenyl)-N-phenyl-2-propenamide
Synonym: Tyrphostin AG-494
MF: C₁₆H₁₂N₂O₃
FW: 280.3
Purity: ≥98%
UV/Vis.: λ_{max}: 370 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-494 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-494 in the solvent of choice, which should be purged with an inert gas. AG-494 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AG-494 in these solvents is approximately 1, 20, and 30 mg/ml, respectively.

AG-494 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AG-494 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AG-494 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Protein tyrosine kinase (PTK) inhibitors are potential antiproliferative agents for diseases caused by the hyperactivity of PTKs. Tyrphostins are a class of antiproliferative compounds that act as PTK blockers. PTK inhibitors that preferentially inhibit the epidermal growth factor (EGF) receptor kinase block EGF-dependent cell proliferation.¹ AG-494 is an inhibitor of EGF receptor kinase with an IC₅₀ value of 1 μM in HT-22 cells.²

References

1. Gazit, A., Yaish, P., Gilon, C., *et al.* Tyrphostins I: Synthesis and biological activity of protein tyrosine kinase inhibitors. *J. Med. Chem.* **32**(10), 2344-2352 (1989).
2. Sagara, Y., Ishige, K., Tsai, C., *et al.* Tyrphostins protect neuronal cells from oxidative stress. *J. Biol. Chem.* **277**(39), 36204-36215 (2002).

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, 05/03/2024

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