

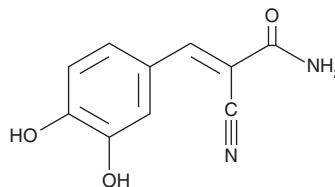
PRODUCT INFORMATION



AG-99

Item No. 10010313

CAS Registry No.: 122520-85-8
Formal Name: 2-cyano-3-(3,4-dihydroxyphenyl)-2E-propenamide
Synonyms: Tyrphostin 46, Tyrphostin AG-99
MF: C₁₀H₈N₂O₃
FW: 204.2
Purity: ≥98%
UV/Vis.: λ_{max}: 254, 364 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-99 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-99 in the solvent of choice, which should be purged with an inert gas. AG-99 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AG-99 in these solvents is approximately 20 mg/ml.

AG-99 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AG-99 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AG-99 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

Protein tyrosine kinase (PTK) inhibitors are potential antiproliferative agents for diseases caused by the hyperactivity of PTKs. Tyrphostins are a class of antiproliferative compounds which act as PTK blockers. PTK inhibitors which preferentially inhibit the epidermal growth factor (EGF) receptor kinase block EGF-dependent cell proliferation.¹ AG-99 is an inhibitor of EGF receptor kinase with an IC₅₀ value of 10 μM in the human epidermoid carcinoma cell line A431.¹

Reference

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).

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, 01/26/2024

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