

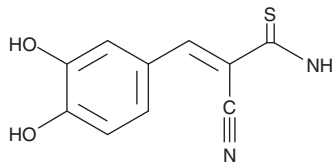
# PRODUCT INFORMATION



## AG-213

Item No. 10010314

**CAS Registry No.:** 122520-86-9  
**Formal Name:** 2-cyano-3-(3,4-dihydroxyphenyl)-2E-propenethioamide  
**Synonyms:** Tyrphostin AG-213, Tyrphostin 47  
**MF:** C<sub>10</sub>H<sub>8</sub>N<sub>2</sub>O<sub>2</sub>S  
**FW:** 220.2  
**Purity:** ≥97%  
**UV/Vis.:** λ<sub>max</sub>: 249, 323, 392 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-213 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-213 in the solvent of choice, which should be purged with an inert gas. AG-213 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AG-213 in these solvents is approximately 10, 20, and 15 mg/ml, respectively.

AG-213 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AG-213 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AG-213 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 that act as PTK blockers. PTK inhibitors which preferentially inhibit the epidermal growth factor (EGF) receptor kinase block EGF-dependent cell proliferation.<sup>1</sup> AG-213 is an inhibitor of EGF receptor kinase with an IC<sub>50</sub> value of 2.4 μM in the human epidermoid carcinoma cell line A431.<sup>1</sup>

### 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, 05/09/2024

#### 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