# **AG-494**

#### **EGFR** inhibitor

AG-494 is a member of the tyrphostin family of tyrosine kinase inhibitors and is a potent inhibitor of EGF receptor autophosphorylation (IC $_{50}$ =1.2  $\mu$ M) and EGF-dependent cell growth (IC $_{50}$ =6  $\mu$ M). It selectively inhibits HER1 (EGF receptor) vs. HER1-2 receptor autophosphorylation (HER1: IC $_{50}$ =1.1  $\mu$ M; HER1-2: IC $_{50}$ =45  $\mu$ M). HER1-2 is a chimeric receptor consisting of the external HER1 domain fused to an internal HER2 domain. Blocks CDK2 activation and causescells to arrest at late G1 and during S phase.

10mg

Citations: 5

**View Online »** 

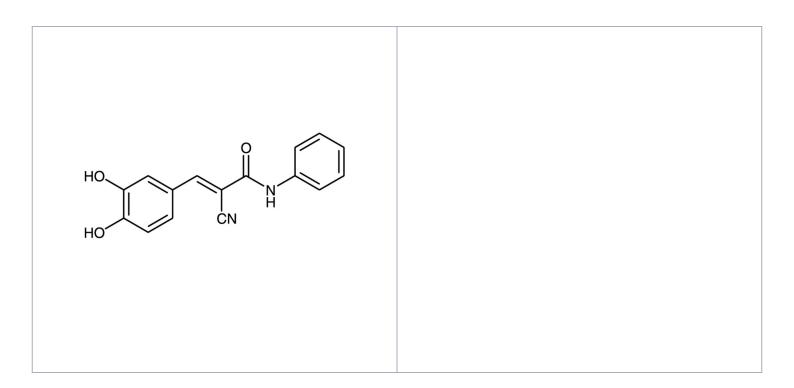
**Ordering Information** 

Order Online »

BML-EI228-0010

Manuals, SDS & CofA

**View Online** »



### **Handling & Storage**

Use/Stability As indicated on product label or CoA when stored as recommended. Product is subject

to oxidation.

**Handling** Store tightly sealed in the dark. Protect from light. Make solutions fresh daily.

Long Term Storage Ambient

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Appearance Yellow solid.

**CAS** 133550-35-3

Couple Target EGFR

Couple Type Inhibitor

Formula  $C_{16}H_{12}N_2O_3$ 

MW 280.3

Purity ≥98% (TLC)

**Solubility** Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).

Last modified: May 29, 2024