

# 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 ( $IC_{50}=1.2\text{ }\mu\text{M}$ ) and EGF-dependent cell growth ( $IC_{50}=6\text{ }\mu\text{M}$ ). It selectively inhibits HER1 (EGF receptor) vs. HER1-2 receptor autophosphorylation (HER1:  $IC_{50}=1.1\text{ }\mu\text{M}$ ; HER1-2:  $IC_{50}=45\text{ }\mu\text{M}$ ). HER1-2 is a chimeric receptor consisting of the external HER1 domain fused to an internal HER2 domain. Blocks CDK2 activation and causes cells to arrest at late G1 and during S phase.

Citations: 5

[View Online »](#)

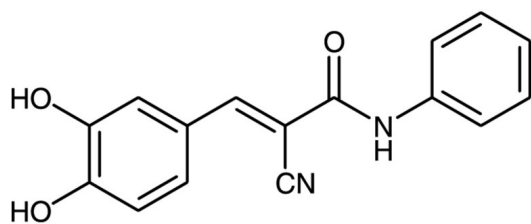
## Ordering Information

[Order Online »](#)

BML-EI228-0010	10mg
----------------	------

## 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).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)