



AG 825

2-Hydroxy-3-methoxy-5-(benzothiazolylthiomethyl)-benzylidenecyanoacetamide

PRODUCT ANALYSIS SHEET

Catalog Number:	PHZ1214
Lot Number:	See product label
Quantity:	2 mg
Appearance:	Yellow solid
Molecular Formula:	C ₁₉ H ₁₅ N ₃ O ₃ S ₂
Molecular Weight:	397.5
Purity:	≥95%, as determined by HPLC
Summary:	AG 825, known alternatively as Tyrphostin AG 825, is a potent, selective inhibitor of ErbB-2 (HER2) autophosphorylation. AG 825 is less effective at inhibiting epidermal growth factor (EGF) receptor autophosphorylation. This compound induces apoptosis with androgen-independent prostate cancer cells, and enhances chemosensitivity with non-small cell lung cancer and ovarian cancer cells.
Biological Activity:	ErbB-2: IC ₅₀ = 0.35 μM Epidermal Growth Factor Receptor: IC ₅₀ = 19 μM
Solubility:	Soluble in DMSO at a concentration of 10 mg/mL.
Sterility:	This product is not sterile.
Storage:	Store, as supplied, at -20°C, protected from light. Upon solubilization, apportion into working aliquots and store at -20°C. Avoid repeated freeze/thaw cycles. Solutions are stable at -20°C for up to two months.
Expiration Date:	Expires one year from the date of receipt when stored as instructed.
Related Products:	ErbB-2 [pY1139] antibody, Cat. # 44-902 ErbB-2 antibody, Cat. # AHO0911 ErbB-2 [pY1248] antibody, Cat. # 44-900 ErbB-2 antibody, FITC conjugate Cat. # AHO0918 ErbB-2 antibody, Cat. # AHO1011 ErbB-2 antibody, biotin conjugate Cat. # AHO0919

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Manufactured under ISO 13485 Quality Standard

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI PHZ1214

(Rev 1.1) (DCC-08-1232)

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.



References:

Osherov, N., et al. (1993) Selective inhibition of the epidermal growth factor and HER2/neu receptors by tyrphostins. *J. Biol. Chem.* 268(15):11134-11142.

Levitzki, A. and A. Gazit (1995) Tyrosine kinase inhibition: an approach to drug development. *Science* 267(5205):1782-1788.

Tsai, C.M., et al. (1996) Enhancement of chemosensitivity by tyrphostin AG825 in high-p185(neu) expressing non-small cell lung cancer cells. *Cancer Res.* 56(5):1068-1074.

Murillo, H., et al. (2001) Tyrphostin AG825 triggers p38 mitogen-activated protein kinase-dependent apoptosis in androgen-independent prostate cancer cells C4 and C4-2. *Cancer Res.* 61(20):7408-7412.

Grunt, T.W. (2003) Tyrphostins and retinoids cooperate during inhibition of in vitro growth of ovarian cancer cells. *Cancer Lett.* 189(2):147-156.

Caution:

Avoid contact with eyes, skin, and mucous membranes. Wear protective clothing when handling this product. Not for human use.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Manufactured under ISO 13485 Quality Standard

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI PHZ1214

(Rev 1.1) (DCC-08-1232)

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

 **invitrogen™**