



SU6656

PRODUCT ANALYSIS SHEET

Catalog Number:	PHZ1203	
Lot Number:	See product label	
Quantity:	1 mg	
Appearance:	Orange solid. Packaged under an inert gas.	
Molecular Formula:	$C_{19}H_{21}N_3O_3S$	
Molecular Weight:	371.5	
Purity:	$\geq 95\%$, as determined by HPLC	
Summary:	SU6656 is a potent inhibitor of Src-family protein tyrosine kinases. Inhibition is competitive with respect to ATP. This compound is useful for elucidating the roles of Src family members in a variety of cell functions, including receptor internalization, motility, apoptosis, differentiation, oncogenesis, cell cycle control, and axonal growth. This compound also has utility in examining the role of Src family members in cellular responses to hyperosmotic stress.	
Biological Activity:	Src: IC_{50} = 280 nM Fyn: IC_{50} = 170 nM Yes: IC_{50} = 20 nM Lyn: IC_{50} = 130 nM	Lck: IC_{50} = 6.88 μ M PDGFR β receptor tyrosine kinase: IC_{50} > 10 μ M IGF-1R tyrosine kinase: IC_{50} > 20 μ M
Solubility:	Soluble in DMSO at a concentration of 25 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 three months.	
Expiration Date:	Expires one year from date of receipt when stored as instructed.	
Related Products:	Src pan antibody, Cat. # 44-655G Src pan antibody, Cat. # 44-656G Fyn pan antibody, Cat. # AHO0652 Fyn pan antibody, Cat. # AHO0482	Src [pY418] antibody, Cat. # 44-660G Src [pY529] antibody, Cat. # 44-662G Src family negative regulatory site (phosphotyrosine) antibody, Cat. # 44-912
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 PHZ1203

(Rev 2.0) 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:

- Bowman, T., et al. (2001) Stat3-mediated Myc expression is required for Src transformation and PDGF-induced mitogenesis. *Proc. Nat'l. Acad. Sci. USA* 98(13):7319-7324.
- Blake, R.A., et al. (2000) SU6656, a selective src family kinase inhibitor, used to probe growth factor signaling. *Mol. Cell Biol.* 20(23):9018-9027.
- Sorkina, T., et al. (2002) Effect of tyrosine kinase inhibitors on clathrin-coated pit recruitment and internalization of epidermal growth factor receptor. *J. Biol. Chem.* 277(30):27433-27441.
- Sanguinetti, A.R., et al. (2003) Fyn is required for oxidative- and hyperosmotic-stress-induced tyrosine phosphorylation of caveolin-1. *Biochem. J.* 376(Pt 1):159-168.
- Reinehr, R., et al. (2004) The Src family kinase Yes triggers hyperosmotic activation of the epidermal growth factor receptor and CD95. *J. Biol. Chem.* 279(23):23977-23987.
- Paveliev, M., et al. (2004) GDNF family ligands activate multiple events during axonal growth in mature sensory neurons. *Mol. Cell Neurosci.* 25(3):453-459.
- Contri, A., et al. (2005) Chronic lymphocytic leukemia B cells contain anomalous Lyn tyrosine kinase, a putative contribution to defective apoptosis. *J. Clin. Invest.* 115(2):369-378.
- Lannutti, B.J., et al. (2005) Induction of polyploidization in leukemic cell lines and primary bone marrow by Src kinase inhibitor SU6656. *Blood* Jan 27; [Epub ahead of print].

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 PHZ1203

(Rev 2.0) 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™**