



PP1

4-amino-5-(4-methylphenyl)-7-(t-butyl)pyrazolo[3,4-d]pyrimidine

PRODUCT ANALYSIS SHEET

| | |
|-----------------------------|--|
| Catalog Number: | PHZ1213 |
| Lot Number: | See product label |
| Quantity: | 1.0 mg |
| Appearance: | Tan solid |
| Molecular Formula: | C ₁₆ H ₁₉ N ₅ |
| Molecular Weight: | 281.4 |
| Purity: | 98%, as assessed by TLC |
| Summary: | PP1 is a highly potent and selective inhibitor of Src family tyrosine kinases. Inhibition of T-cell phosphorylation in response to anti-CD3 activation occurs at 10 μM. |
| Biological Activity: | Lck: IC ₅₀ = 5 nM Lyn: IC ₅₀ = 6 nM Src: IC ₅₀ = 170 nM |
| Solubility: | Soluble in DMSO at a concentration of 25 mg/mL. |
| Sterility: | This product is not sterile. |
| Storage: | Store, as supplied, at -20°C. 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: | Lck [pS ¹⁵⁸] antibody, Cat. # 44-842 Src [pY418] antibody, Cat. # 44-660G Lck [pY ¹⁹²] antibody, Cat. # 44-844G Src Family Neg. Reg. [pY] Site antibody, Cat. # 44-912 Lck [pY ⁵⁰⁵] antibody, Cat. # 44-850G Hck [pYpS ^{209/211}] antibody, Cat. # 44-910 |
| References: | Hanke, J.H., et al. (1996) Discovery of a novel, potent, and Src family-selective tyrosine kinase inhibitor. Study of Lck- and FynT-dependent T cell activation. <i>J. Biol. Chem.</i> 271:695-701. DeSouza, N., et al. (2002) Protein kinase A and two phosphatases are components of the inositol 1,4,5-trisphosphate receptor macromolecular signaling complex. <i>J. Biol. Chem.</i> 277(42):39397-39400. Shah, O.J., et al. (2002) The Src-family tyrosine kinase inhibitor PP1 interferes with the activation of ribosomal protein S6 kinases. <i>Biochem. J.</i> 366(Pt 1):57-62. Cohen, P.T. (2002) Protein phosphatase 1: targeted in many directions. <i>J. Cell. Sci.</i> 115(Pt 2):241-256. |
| 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 PHZ1213

(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.

