


PP2 4-amino-5-(4-chlorophenyl)-7-(t-butyl)pyrazolo [3,4-d]pyrimidine

Publication Number MAN0003710

Rev. 2.00








Catalog Number:	PHZ1223
Quantity:	1.0 mg
Lot Number:	See product label.
Appearance:	Off-white solid
Molecular Formula:	C ₁₅ H ₁₆ ClN ₅
Molecular Weight:	301.8
Purity:	98%, as assessed by TLC.
Summary:	PP2 is a potent and selective inhibitor of the Src family tyrosine kinases. PP2 is a useful tool for probing the involvement of Src family tyrosine kinases in cellular signaling.
Biological Activity:	Lck: IC ₅₀ = 4 nM Hck: IC ₅₀ = 5 nM Fyn: IC ₅₀ = 5 nM EGFR: IC ₅₀ = 480 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.
	Avoid contact with eyes, skin, and mucous membranes. Wear protective clothing when handling this product. Not for human use.





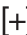

Related Products:

Product	Catalog no.
EGFR Sample Pack	44-799G
Lck [pS ¹⁵⁸] antibody	44-842
Lck [pY ¹⁹²] antibody	44-844G
Lck [pY ⁵⁰⁵] antibody	44-850G
Src [pY418] antibody	44-660G
Src Family Neg. Reg. [pY] Site antibody	44-912
Hck [pYpS ^{209/211}] antibody	44-910
FAK [pY ⁸⁶¹] antibody	44-626G

Explanation of Symbols

The symbols present on the product label are explained below:

Symbol	Description
	Catalog Number
	Research Use Only
	Use by
	Manufacturer
	Without, does not contain
	Protect from light
	Directs the user to consult instructions for use (IFU), accompanying the product.

Symbol	Description
	Batch code
	In vitro diagnostic medical device
	Temperature limitation
	European Community authorized representative
	With, contains
	Consult accompanying documents

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

For Research Use Only. Caution: Not for human or animal therapeutic or diagnostic use.

Manufacturing site: 7335 Executive Way | Frederick, MD 21704 | Toll Free in USA 800.955.6288

© 2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

Revision Date 13 May 2011

