



**p38 MAPK Inhibitor**  
2-(4-Chlorophenyl)-4-(4-fluorophenyl)-5-pyridin-4-yl-1,2-dihydropyrazol-3-one

**PRODUCT ANALYSIS SHEET**

<b>Catalog Number:</b>	PHZ1112
<b>Lot Number:</b>	See product label
<b>Quantity:</b>	0.5 mg
<b>Appearance:</b>	Orange-brown solid. Packaged under an inert gas.
<b>Molecular Formula:</b>	C <sub>20</sub> H <sub>13</sub> ClFN <sub>3</sub> O
<b>Molecular Weight:</b>	365.8
<b>Purity:</b>	≥95%, as determined by HPLC
<b>Summary:</b>	This compound is a potent inhibitor of p38 MAPK. Inhibition of PKC, PKG, and MLCK is negligible. When applied to cultured neurons, this compound is observed to inhibit axon branching.
<b>Biological Activity:</b>	p38 MAPK: IC <sub>50</sub> = 35 nM
<b>Solubility:</b>	Soluble in DMSO at a concentration of 25 mg/mL.
<b>Sterility:</b>	This product is not sterile.
<b>Storage:</b>	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.
<b>Expiration Date:</b>	Expires one year from date of receipt when stored as instructed.
<b>Related Products:</b>	p38 MAPK [pTpY180/182] antibody, Cat. # 44-684G p38 MAPK antibody, Cat. # AHO0782 ERK1&2 [pTpY185/187] antibody, Cat. # 44-680G JNK1&2 [pTpY183/185] antibody, Cat. # 44-682G
<b>Reference:</b>	de Laszlo, S.E., et al. (1998) Pyrroles and other heterocycles as inhibitors of p38 kinase. Bioorg. Med. Chem. Lett. 8(19):2689-2894.
<b>Caution:</b>	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](http://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](mailto:techsupport@invitrogen.com)

PI PHZ1112

(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](http://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™