

Catalog Number:	PHZ1271												
Lot Number:	See product label												
Quantity:	100 µg												
Appearance:	Lyophilized solid												
Molecular Formula:	C ₂₈ H ₂₆ N ₄ O ₃												
Molecular Weight:	466.5												
Purity:	99%, as assessed by TLC												
Summary:	Staurosporine inhibits a variety of kinases including PKA, PKG, MLCK, CaMK, tyrosine kinases, and phosphorylase kinase. This compound displays selectivity toward the PKC isoforms, inhibiting PKCα, PKCβ, PKCγ, PKCδ, and PKCε, but not PKCζ, and induces PKC translocation. Staurosporine also inhibits topoisomerase II directly by interaction with the ATP binding site, augments PMA-induced ornithine decarboxylase, and activates a bcl-2-regulated apoptosis pathway.												
Biological Activity:	<table> <tr> <td>CaMK: IC₅₀ = 20 nM</td><td>PKC: K_i = 0.7 nM</td></tr> <tr> <td>MLCK: K_i = 1.3 nM</td><td>PKCα: IC₅₀ = 58 nM</td></tr> <tr> <td>phosphorylase kinase IC₅₀ = 0.5 nM</td><td>PKCβ: IC₅₀ = 65 nM</td></tr> <tr> <td>PKA: K_i = 7.0 nM</td><td>PKCγ: IC₅₀ = 49 nM</td></tr> <tr> <td>PKG: K_i = 8.5 nM</td><td>PKCδ: IC₅₀ = 325 nM</td></tr> <tr> <td>tyrosine kinases: IC₅₀ = 70 nM</td><td>PKCε: IC₅₀ = 160 nM</td></tr> </table>	CaMK: IC ₅₀ = 20 nM	PKC: K _i = 0.7 nM	MLCK: K _i = 1.3 nM	PKCα: IC ₅₀ = 58 nM	phosphorylase kinase IC ₅₀ = 0.5 nM	PKCβ: IC ₅₀ = 65 nM	PKA: K _i = 7.0 nM	PKCγ: IC ₅₀ = 49 nM	PKG: K _i = 8.5 nM	PKCδ: IC ₅₀ = 325 nM	tyrosine kinases: IC ₅₀ = 70 nM	PKCε: IC ₅₀ = 160 nM
CaMK: IC ₅₀ = 20 nM	PKC: K _i = 0.7 nM												
MLCK: K _i = 1.3 nM	PKCα: IC ₅₀ = 58 nM												
phosphorylase kinase IC ₅₀ = 0.5 nM	PKCβ: IC ₅₀ = 65 nM												
PKA: K _i = 7.0 nM	PKCγ: IC ₅₀ = 49 nM												
PKG: K _i = 8.5 nM	PKCδ: IC ₅₀ = 325 nM												
tyrosine kinases: IC ₅₀ = 70 nM	PKCε: IC ₅₀ = 160 nM												
Solubility:	Soluble in DMSO at a concentration of 25 mg/mL and DMF at a concentration of 25 mg/mL.												
Storage:	Store, as supplied, at 2-8°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:	<table> <tr> <td>ERK1&2 [pTpY^{185/187}] antibody, Cat. # 44-680G</td><td>Tau [pS²¹⁴] antibody, Cat. # 44-742G</td></tr> <tr> <td>Akt/PKB [pS⁴⁷³] antibody, Cat. # 44-622G</td><td>Tau [pS²⁶²] antibody, Cat. # 44-750G</td></tr> <tr> <td>PKCδ [pY³¹¹] (mouse) antibody, Cat. # 44-950</td><td>Tau [pS³⁵⁶] antibody, Cat. # 44-751G</td></tr> <tr> <td>Tau [pT²¹²] antibody, Cat. # 44-740G</td><td>Tau [pS⁴⁰⁹] antibody, Cat. # 44-760G</td></tr> </table>	ERK1&2 [pTpY ^{185/187}] antibody, Cat. # 44-680G	Tau [pS ²¹⁴] antibody, Cat. # 44-742G	Akt/PKB [pS ⁴⁷³] antibody, Cat. # 44-622G	Tau [pS ²⁶²] antibody, Cat. # 44-750G	PKCδ [pY ³¹¹] (mouse) antibody, Cat. # 44-950	Tau [pS ³⁵⁶] antibody, Cat. # 44-751G	Tau [pT ²¹²] antibody, Cat. # 44-740G	Tau [pS ⁴⁰⁹] antibody, Cat. # 44-760G				
ERK1&2 [pTpY ^{185/187}] antibody, Cat. # 44-680G	Tau [pS ²¹⁴] antibody, Cat. # 44-742G												
Akt/PKB [pS ⁴⁷³] antibody, Cat. # 44-622G	Tau [pS ²⁶²] antibody, Cat. # 44-750G												
PKCδ [pY ³¹¹] (mouse) antibody, Cat. # 44-950	Tau [pS ³⁵⁶] antibody, Cat. # 44-751G												
Tau [pT ²¹²] antibody, Cat. # 44-740G	Tau [pS ⁴⁰⁹] antibody, Cat. # 44-760G												
Caution:	Avoid contact with eyes, skin, and mucous membranes. Wear protective clothing when handling this product. Not for human use.												
References:	<p>Seynaeve, C.M., et al. (1994) Differential inhibition of protein kinase C isozymes by UCN-01, a staurosporine analogue. <i>Mol. Pharmacol.</i> 45(6):1207-1214.</p> <p>Wang, X., et al. (1996) Cleavage of sterol regulatory element binding proteins (SREBPs) by CPP32 during apoptosis. <i>EMBO J.</i> 15(5):1012-1020.</p> <p>Spyridopoulos, I., et al. (2002) Divergence of angiogenic and vascular permeability signaling by VEGF: inhibition of protein kinase C suppresses VEGF-induced angiogenesis, but promotes VEGF-induced, NO-dependent vascular permeability. <i>Arterioscler. Thromb. Vasc. Biol.</i> 22(6):901-906.</p> <p>Wan, X., et al. (2002) PTEN augments staurosporine-induced apoptosis in PTEN-null Ishikawa cells by downregulating PI3K/Akt signaling pathway. <i>Cell Death Differ.</i> 9(4):414-420.</p> <p>Swannie, H.C. and S.B. Kaye (2002) Protein kinase C inhibitors. <i>Curr. Oncol. Rep.</i> 4(1):37-46.</p>												

For research use only. CAUTION: Not intended for human or animal therapeutic or diagnostic use.

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 PHZ1271

(Rev 4.0) DCC-09-1799

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.