



KT5720

PRODUCT ANALYSIS SHEET

Catalog Number:	PHZ1141
Lot Number:	See product label
Quantity:	0.1 mg
Appearance:	White solid
Molecular Formula:	$C_{32}H_{31}N_3O_5$
Molecular Weight:	537.6
Purity:	≥95%, as determined by HPLC
Summary:	KT5720 is a potent, cell permeable inhibitor of PKA, and is a valuable tool in studies delineating cAMP-dependent signaling pathways, including activation of the transcription factor CREB. Inhibition of PKC, PKG, and MLCK is negligible. KT5720 is prepared by a chemical modification of K-252a. This compound has utility in neuroscience studies, and has been used to examine the role PKA in CNTF-induced neuronal survival, BDNF-induced neuronal plasticity, and plasminogen and ganglioside-induced activation of microglia. This compound has also been used to examine the role of PKA in modulating the activity of ion channels and synaptic transmission.
Biological Activity:	PKA: $K_i = 56$ nM
Solubility:	Soluble in DMSO at a concentration of 1 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 six months.
Expiration Date:	Expires one year from date of receipt when stored as instructed.
Related Products:	PKA α/β [pT197] antibody, Cat. # 44-988 PKA β [pS338] antibody, Cat. # 44-992G PKA II β [pS114] antibody, Cat. # 44-998G
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 PHZ1141

(Rev 1.1) (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™

References:

- Kase, H., et al. (1987) K-252 compounds, novel and potent inhibitors of protein kinase C and cyclic nucleotide-dependent protein kinases. *Biochem. Biophys. Res. Commun.* 142(2):436-440.
- Gadbois, D.M., et al. (1992) Multiple kinase arrest points in the G1 phase of nontransformed mammalian cells are absent in transformed cells. *Proc. Nat'l. Acad. Sci.* 89(18):8626-8630.
- Cabell, L. and G. Audesirk (1993) Effects of selective inhibition of protein kinase C, cyclic AMP-dependent protein kinase, and Ca(2+)-calmodulin-dependent protein kinase on neurite development in cultured rat hippocampal neurons. *Int. J. Dev. Neurosci.* 11(3):357-368.
- Simpson, C.S. and B.J. Morris (1995) Induction of c-fos and zif/268 gene expression in rat striatal neurons, following stimulation of D1-like dopamine receptors, involves protein kinase A and protein kinase C. *Neuroscience* 68(1):97-106.
- Ellershaw, D.C., et al. (2002) Modulation of volume-sensitive chloride current by noradrenaline in rabbit portal vein myocytes. *J. Physiol.* 542(Pt 2):537-547.
- Kumar, A., et al. (2002) Distinct signaling pathways are activated in response to mechanical stress applied axially and transversely to skeletal muscle fibers. *J. Biol. Chem.* 277(48):46493-46503.
- Blanquet, P.R., et al. (2003) A calcium/calmodulin kinase pathway connects brain-derived neurotrophic factor to the cyclic AMP-responsive transcription factor in the rat hippocampus. *Neuroscience* 118(2):477-490.
- Park, K., et al. (2004) Cellular mechanisms associated with spontaneous and ciliary neurotrophic factor-cAMP-induced survival and axonal regeneration of adult retinal ganglion cells. *Neuroscience* 24(48):10806-10815.
- Min, K.J., et al. (2004) Protein kinase A mediates microglial activation induced by plasminogen and gangliosides. *Exp. Mol. Med.* 36(5):461-467.

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 PHZ1141

(Rev 1.1) (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™