



ERK Activation Inhibitor Peptide I, Cell Permeable

PRODUCT ANALYSIS SHEET

Catalog Number:	PHZ1103
Lot Number:	See product label
Appearance:	Lyophilized solid. Packaged under an inert gas.
Form:	Trifluoroacetate salt
Amino Acid Sequence:	Stearoyl-Met-Pro-Lys-Lys-Lys-Pro-Thr-Pro-Ile-Gln-Leu-Asn-Pro-NH ₂
Molecular Weight:	1757.3
Purity:	≥95%, as determined by HPLC
Summary:	ERK Activation Inhibitor Peptide I is a steared, cell permeable peptide that corresponds to the N terminus of MEK1 (MAP kinase kinase 1). This peptide specifically inhibits the activation of ERK2 by blocking its binding to MEK. ERK Activation Inhibitor Peptide I is observed to inhibit ERK activation by PMA or NGF, and to inhibit activation of the transcription factor Elk-1. The activation of JNK and p38 MAPK is not affected by this peptide.
Biological Activity:	ERK2: IC ₅₀ = 2.5 μM
Solubility:	Soluble in water at a concentration of 1 mg/mL or DMSO at a concentration of 5 mg/mL.
Sterility:	This product is not sterile.
Storage:	Store, as supplied, at -20°C, preferably desiccated. Protect 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.
Expiration Date:	Expires one year from date of receipt when stored as instructed.
Related Products:	ERK1/2 [pTpY185/187] antibody, Cat. # 44-680G p38 MAPK [pTpY180/182] antibody, Cat. # 44-684G JNK1/2 [pTpY183/185] antibody, Cat. # 44-682G
References:	Kelemen, B.R., et al. (2002) Selective in vivo inhibition of mitogen-activated protein kinase activation using cell-permeable peptides. J. Biol. Chem. 277(10):8741-8748.
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 PHZ1103

(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.

