

PAKtide

Synthetic peptide substrate

Catalog # P08-58

Lot # L242-3

Product Description

The 10 amino acids of PAKtide peptide (RRRLSFAEPG) contain a serine/threonine protein kinase phosphorylation site in a common seven-residue epitope (1, 2).

Molecular Weight

The theoretical molecular weight is 1188.4

Purity

The purity was determined to be more than 96.9% by HPLC analysis.

Formulation

1mg of peptide supplied as a lyophilized powder.

Storage and Stability

Store product at -20°C. For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Reconstitution Protocol

Dilute peptide in 20mM Tris-HCl, pH 7.5 solution to a final concentration of 1mg/ml.

References

1. Heike ROSS et al: A non-radioactive method for the assay of many serine/threonine-specific protein kinases. *Biochem. J.* (2002) 366, 977±981
2. Hao Wu et al: The Mechanism of p21-activated Kinase 2 Autoactivation. *J. Biol. Chem.*, Vol. 278, Issue 43, 41768-41778, October 24, 2003

Related Kinases

PAKtide can be utilized as a substrate for the following active protein kinases.

Product Name

PAK1, active

Catalog Number

P02-18G

PAKtide

Synthetic peptide substrate

Catalog Number P08-58

Quantity 1 mg

Specific Lot Number L242-3

Purity 96.9%

Format 1 mg lyophilized powder

Stability 1yr At -20°C from date of shipment

Storage & Shipping Store product at -20°C. For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped at ambient temperature.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.