

Histone H1 Peptide (152-159)

Synthetic peptide substrate derived from human Histone H1

Catalog # H10-58

Lot # G1309-6

Product Description

The Histone H1 Peptide (152-159) sequence (GGGPATP-KKAKKL-COOH) is derived from human histone H1 between amino acids 152-159 and is suitable for use as the substrate for CDK family kinase assays, such as CDK1, CDK2, CDK5 and CDK6.

Molecular Weight

The theoretical molecular weight is 1252.53.

Purity

The purity was determined to be 95.19% 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 suitable concentration.

Histone H1 Peptide (152-159)

Synthetic peptide substrate derived from human Histone H1

Catalog #	H10-58
Quantity	1 mg
Lot #	G1309-6
Purity	95.19%
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.