

Histone H4 Peptide (1-21)

Synthetic peptide substrate derived from human Histone H4

Catalog # H13-58

Lot # I292-8

Product Description

The Histone H4 Peptide (1-21) sequence (SGRGKGGKGL-GKGGAKRHRKVGGKKC) is derived from human histone H4 (1-21) and is suitable for use as the substrate for histone methyltransferase (at R3) and acetyltransferase (at K5, K8, K12 and K16) family assays.

Molecular Weight

The theoretical molecular weight is 2565.06

Purity

The purity was determined to be 97.27% by HPLC analysis.

Formulation

0.5mg 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.

Related Kinases

Histone H4 Peptide (1-21) can be utilized as a substrate for histone methyltransferase (at R3) and acetyltransferase (at K5, K8, K12 and K16) family assays. Please see the available related enzyme products at (<http://www.signalchem.com/products.php>).

Histone H4 Peptide (1-21)

Synthetic peptide substrate derived from human Histone H4

Catalog Number	H13-58
Quantity	0.5 mg
Specific Lot Number	I292-8
Purity	97.27%
Format	0.5 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.