

## Acetylated Histone H4 Peptide I

Synthetic peptide substrate derived from human Histone H4

**Catalog # H13-358**

Lot # I292-9

### Product Description

The Acetylated Histone H4 (K5, 8, 12, 16) Peptide sequence (SGRG[Ac-K]GG[Ac-K]GLG[Ac-K]GGA[Ac-K]RHRKVGGKKC) is derived from human histone H4 (1-21) and is suitable for use as the assay control of histone (de-)methylation and (de-)acetylation.

### Molecular Weight

The theoretical molecular weight is 2733.06

### Purity

The purity was determined to be 95.35% 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

Acetylated Histone H4 (K5, 8, 12, 16) Peptide can be utilized as the assay control of histone (de-)methylation and (de-)acetylation. Please see the available related products at <http://www.signalchem.com/products.php>.

## Acetylated Histone H4 Peptide I

Synthetic peptide substrate derived from human Histone H4

Catalog Number	H13-358
Quantity	0.5 mg
Specific Lot Number	I292-9
Purity	95.35%
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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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