

PRODUCT INFORMATION



Histone H4 Peptide Substrate (1-21)

Item No. 10854

Overview and Properties

Peptide Sequence: SGRGKGGKGLGKGGAKRHRKV
Synonym: H4 Peptide
Storage: -20°C (as supplied)
Stability: As supplied, 1 year from the QC date provided on the Certificate of Analysis, when stored properly
Supplied as: A lyophilized peptide
Solubility: At least 20 mg/ml with water or aqueous buffer

Description

Histone H4 undergoes many modifications including acetylation, methylation, and phosphorylation that are important for regulation of gene transcription.¹ Histone H4 peptide can be used as a target substrate for several of the histone modifying enzymes including lysine methyltransferases, arginine methyltransferases, acetyltransferases, and others.

Reference

1. Bhaumik, S.R., Smith, E., and Shilatifard, A. Covalent modifications of histones during development and disease pathogenesis. *Nat. Struct. Mol. Biol.* **14(11)**, 1008-1016 (2007).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/05/2016

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM