

## Histone H1 Protein

Native proteins purified from bovine thymus

**Catalog # H10-54N**

Lot # M119-3

### Product Description

Native histone H1 was purified from bovine thymus tissues. This product contains lysine-rich fraction and is routinely used as a substrate of CDK family.

### Gene Aliases

None

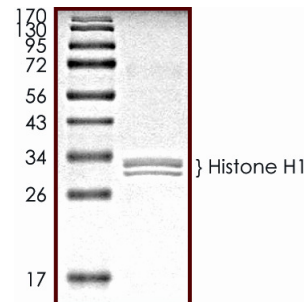
### Formulation

Native protein stored in 50mM Tris-HCl, pH 7.5, and 150mM NaCl.

### Storage and Stability

Store product at  $-20^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

### Purity



The purity of Histone H1 was determined to be **>95%** by densitometry. The major protein band was approx. MW **31~33kDa**.

## Histone H1 Protein

Native proteins purified from bovine thymus

Catalog # H10-54N

Lot # M119-3

Purity >95%

Concentration 1  $\mu\text{g}/\mu\text{l}$

Stability 2yr at  $-20^{\circ}\text{C}$  from date of shipment

Storage & Shipping Store product at  $-20^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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.**