

## Histone H3 Protein

Native proteins purified from bovine thymus

Catalog # H12-54N

Lot # X695-1

### Product Description

Native histone H3 was purified from bovine thymus tissues. This product contains lysine-rich fraction and is routinely used as a substrate of SLK kinase.

### Gene Aliases

None

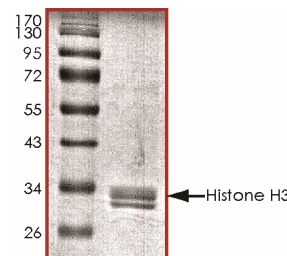
### Formulation

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 H3 was determined to be **>95%** by densitometry.

The major protein bands were approx. MW **30~34kDa**.

## Histone H3 Protein

Native proteins purified from bovine thymus

Catalog # H12-54N

Lot # X695-1

Purity >95%

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

Stability 1yr 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.