

Histone H1 Protein

Native proteins purified from bovine thymus

Catalog # H10-54N

Lot # R182-3

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

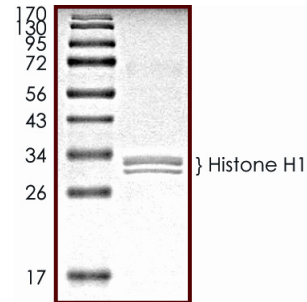
Formulation

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

Storage and Stability

Store product at -20°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 # R182-3

Purity >95%

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

Stability 2yr at -20°C from date of shipment

Storage & Shipping Store product at -20°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.