

## HeLa Core Histones

---

**Catalog No:** 53501

**Contents:** 24 µl of HeLa Core Histones at 1.5 µg/µl

**Composition:**

HeLa Core Histones were purified from HeLa chromatin by hydroxylapatite column chromatography and are formulated in a storage buffer containing 10 mM HEPES pH 7.9, 1 mM EDTA, 10 mM KCl, 10% glycerol, 0.5 mM PMSF, and 1 mM DTT.

**Quality Control:**

HeLa Core Histones are tested in the Chromatin Assembly Kit, Catalog No. 53500.

**Storage:**

-80°C

**Guarantee:**

This product is guaranteed stable for six months when stored properly.

