

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.

