

## HeLa nuclear extract

---

**Catalog No:** 36010

**Contents:** 2 x 100 µg of HeLa nuclear extract at 2.5 µg/µl

**Composition:**

Nuclear extracts are supplied in lysis buffer (20mM Hepes (pH 7.5), 350 mM NaCl, 20% glycerol, 1% Igepal, 1mM MgCl<sub>2</sub>, 0.1mM EGTA, 5mM DTT and protease inhibitors.

**Quality Control:**

Extracts have been quality control tested by Western blot and TransAM Transcription Factor ELISAs.

**Storage:**

To ensure stability, extracts should be stored at -80°C.

We recommend aliquoting the extracts into single-use fractions and then storing them at -80°C. This eliminates repeated freeze/thaw cycles.

**Guarantee:**

This product is guaranteed for 6 months from date of arrival.