

## HeLa nuclear extract (Anisomycin treated)

---

**Catalog No:** 36111

**Contents:** 2 x 100 µg of HeLa nuclear extract (Anisomycin treated) at 2.5 µg/µl

**Composition:**

Nuclear extracts are supplied in dilution buffer (20 mM Hepes pH 7.5, 400 mM NaCl, 20% glycerol, 0.1 mM EDTA, 10 mM NaF, 10 µM Na<sub>2</sub>MoO<sub>4</sub>, 1 mM NaVO<sub>3</sub>, 10mM PNPP, 10 mM β-glycerophosphate, 1 mM DTT and protease inhibitors). Cells were cultured in medium supplemented with 5 µM Anisomycin for 1 hour at 37°C immediately prior to harvesting.

**Quality Control:**

Each lot has been tested for ATF-2 activation by using TransAM™ Kits.

**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.