

## LoVo nuclear extract

---

**Catalog No:** 36097

**Contents:** 200 µg of LoVo nuclear extract at 7.19 µg/µl

**Composition:**

Nuclear extracts are supplied in dilution buffer (20mM Hepes (pH 7.9), 100 mM KCl, 1mM MgCl<sub>2</sub>, 20% glycerol, 0.5 mM PMSF and 0.5 mM DTT).

**Quality Control:**

Extracts have been quality control tested by Western blot and the Electrophoretic Mobility Shift Assay (EMSA).

**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.