

## Nb2 nuclear extract (prolactin stimulated)

---

**Catalog No:** 36106

**Contents:** 2 x 100 µg of Nb2 nuclear extract (prolactin stimulated) at 2.5 µg/µl

**Composition:**

The Nb2 nuclear extract derived from rat lymphocytes have been collected in Lysis Buffer after a 20 minute stimulation with 25 ng/ml of prolactin at 37°C immediately prior to harvesting. Lysis Buffer consists of 20 mM Hepes pH 7.9, 100 mM KCl, 20% glycerol, 1 mM MgCl<sub>2</sub>, 0.5 mM DTT and 0.5 mM PMSF. The protein content has been determined by a Bradford-based assay.

**Quality Control:**

Each lot has been tested for STAT activation by using TransAM™ STAT Family 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.