

## Jurkat nuclear extract (PHA stimulated)

---

**Catalog No:** 36108

**Contents:** 2 x 100 µg of Jurkat nuclear extract (PHA stimulated) at 2.5 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). Cells were cultured in medium supplemented with 2 µg/ml PHA (Leucoagglutinin, PHA-L) for 2 days at 37°C immediately prior to harvesting.

**Quality Control:**

Each lot has been tested by WB and for NFATc1 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.