

## THP-1 nuclear extract (PMA stimulated)

---

**Catalog No:** 40700

**Contents:** 2 x 100 µg of THP-1 nuclear extract (PMA stimulated) at 2.50 µg/µl

**Composition:**

The THP-1 nuclear extracts have been collected in Lysis Buffer after a 24-hour incubation with PMA (10 ng/ml). The Lysis Buffer consists of 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. The protein content has been determined by a Bradford-based assay.

**Quality Control:**

Extracts have been quality control tested by Western blot.

**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.