

## COS-7 nuclear extract (IFN $\gamma$ treated)

---

**Catalog No:** 36112

**Contents:** 2 x 100  $\mu$ g of COS-7 nuclear extract (IFN $\gamma$  treated) at 2.5  $\mu$ g/ $\mu$ l

**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  $\mu$ M Na<sub>2</sub>MoO<sub>4</sub>, 1 mM NaVO<sub>3</sub>, 10mM PNPP, 10 mM  $\beta$ -glycerophosphate, 1 mM DTT and protease inhibitors). Cells were treated with 1000 U/ml of IFN $\gamma$  for 15 minutes immediately prior to harvesting.

**Quality Control:**

Each lot has been tested for STAT 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.