

## MCF-7 nuclear extract (H<sub>2</sub>O<sub>2</sub> treated)

---

**Catalog No:** 40800

**Contents:** 2 x 100 µg of MCF-7 nuclear extract (H<sub>2</sub>O<sub>2</sub> treated) at 2.50 µg/µl

### **Composition:**

The MCF-7 nuclear extracts have been collected in Lysis Buffer after 3 hrs incubation with 0.2 mM H<sub>2</sub>O<sub>2</sub>. The Lysis Buffer consists of 20mM Hepes (pH 7.9), 100 mM KCl, 1mM MgCl<sub>2</sub>, 20% glycerol, 0.5 mM PMSF and 0.5 mM DTT. The protein content has been determined by a Bradford-based assay.

### **Quality Control:**

Each lot has been tested for p53 activation by using TransAM™ p53 Kit. The signal intensity for p53 activation in each lot is compared to the signal intensity obtained with extracts from untreated MCF-7 cells (see figure). Once the signals are blanked, the ratio between the signals from extracts of treated cells over untreated cells used at 5 µg/well must be above 4. This ratio may vary depending of basal level of p53 activation in a given cell type.

### **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.