

MCF-7 nuclear extract (H₂O₂ post-treated)

Catalog No: 40810

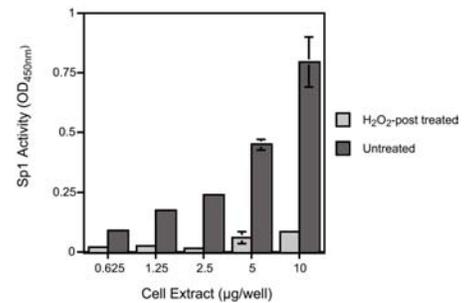
Contents: 2 x 100 µg of MCF-7 nuclear extract (H₂O₂ post treated) at 2.50 µg/µl

Composition:

The MCF-7 nuclear extracts have been collected in Lysis Buffer. Post-collection treatment consists in exposing the nuclear extract to a final concentration of 200 mM H₂O₂ for 20 min on ice. The Lysis Buffer consists of 20mM Hepes (pH 7.9), 100 mM KCl, 1mM MgCl₂, 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 Sp1 activity by using the TransAM™ Sp1 Kit. The signal intensity for Sp1 activity in each lot is compared to the signal intensity obtained with H₂O₂-post treated extracts from MCF-7 cells (see figure). Once the signals are blanked, the ratio between the signals from untreated extracts over treated extracts used at 5 µg/well must be above 4.



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.