

HeLa nuclear extract (TPA stimulated, 2 hr)

Catalog No: 36009

Contents: 2 x 100 µg of HeLa nuclear extract (TPA stimulated, 2 hr) at 2.5 µg/µl

Composition:

Nuclear extracts are supplied in dilution buffer (20mM Hepes (pH 7.9), 100 mM KCl, 1mM MgCl₂, 20% glycerol, 0.5 mM PMSF and 0.5 mM DTT). Cells were cultured in medium supplemented with 50 ng/mL TPA (phorbol, 12-myristate, 13 acetate) for two hours at 37°C immediately prior to harvesting.

Quality Control:

Extracts have been quality control tested by Western blot and TransAM Transcription Factor ELISAs.

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.