

Hep G2 nuclear extract (Acetaldehyde treated)

Catalog No: 36065

Contents: 2 x 100 µg of Hep G2 nuclear extract (Acetaldehyde treated) 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 200 µM Acetaldehyde for 4 hours at 37°C immediately prior to harvesting.

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.