

A-431 nuclear extract (EGF treated)

Catalog No: 36003

Contents: 200 µg of A-431 nuclear extract (EGF treated)

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 10 ng/mL EGF (epidermal growth factor) for 10 minutes 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.

Product Guarantee

This product is guaranteed for 6 months from date of arrival.