

A-431 nuclear extract

Catalog No: 36004

Contents: 200 µg of A-431 nuclear extract

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).

Quality Control

Extracts have been quality control tested by Western blot and the Electrophoretic Mobility Shift Assay (EMSA).

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.