

Mouse Brain nuclear extract

Catalog No: 36053

Contents: 120 µg of Mouse Brain nuclear extract at 5.44 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).

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.

Guarantee:

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