

COS-7 nuclear extract

Catalog No: 36079

Contents: 2 x 100 µg of COS-7 nuclear extract at 2.5 µg/µl

Composition:

Nuclear extracts are supplied in lysis buffer (20mM Hepes (pH 7.5), 400 mM NaCl, 20% glycerol, 0.1 mM EDTA, 10 mM NaF, 10 µM Na₂MoO₄, 1 mM NaVO₃, 10 mM PNPP, 10 mM beta-glycerophosphate, 1 mM DTT and protease inhibitors).

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.