

PC-12 nuclear extract

Catalog No: 36022

Contents: 2 x 100 µg of PC-12 nuclear extract at 2.5 µg/µl

Composition:

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

Quality Control:

Extracts have been quality control tested by Western blot and the TransAM Transcription Factor ELISA Kit.

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.