

PC-12 nuclear extract (Forskolin stimulated)

Catalog No: 40400

Contents: 2 x 100 µg of PC-12 nuclear extract (Forskolin stimulated) at 2.50 µg/µl

Composition:

The PC12 nuclear extracts have been collected in Lysis Buffer after a 30-minute incubation with forskolin (20 µM). The Lysis Buffer consists of 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. The protein content has been determined by a Bradford-based assay.

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.