

WI-38 nuclear extract (Forskolin stimulated)

Catalog No: 40300

Contents: 2 x 100 µg of WI-38 nuclear extract (Forskolin stimulated) at 2.50 µg/µl

Composition:

The WI-38 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, 10 mM PNPP, 10 mM β-glycerophosphate, 1 mM DTT and protease inhibitors. The protein content has been determined by a Bradford-based assay.

Quality Control:

Each lot has been tested for CREB activation by using TransAM™ pCREB Kits. The signal intensity for CREB activation in each lot is compared to the signal intensity obtained with extracts from unstimulated WI-38 VA13 cells (see figure). After the signals are blanked, the ratio of the signals from stimulated cells over unstimulated cells must be above 3. This ratio may vary depending on the basal level of CREB activation in a given cell type.

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.