

WI-38 nuclear extract

Catalog No: 40310

Contents: 2 x 100 µg of WI-38 nuclear extract at 2.5 µg/µl

Composition:

The WI-38 nuclear extracts have been collected in Lysis Buffer. 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:

Each lot has been tested for pCREB activation by using TransAM™ pCREB Kits. The signal intensity for pCREB 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 2 when using TransAM pCREB. This ratio may vary depending on the basal level of pCREB 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.