

WI-38 nuclear extract (TPA + CI stimulated)

Catalog No: 40500

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

Composition:

The WI-38 nuclear extracts have been collected in Lysis Buffer after a 2-hour incubation with TPA (100 ng/ml) and calcium ionophore (2 µg/ml). 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 AP-1 activation by using TransAM™ AP-1 Family Kits. Incubation with wild-type competitor oligos reduces binding by over 80% while incubation with mutant oligos had no effect on AP-1 Family binding to DNA (see figure).

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.