

COS-7 nuclear extract (CoCl₂ treated)

Catalog No: 40600

Contents: 2 x 100 µg of COS-7 nuclear extract (CoCl₂ treated) at 2.50 µg/µl

Composition:

The COS-7 nuclear extracts have been collected in Lysis Buffer after a 16-hour incubation with CoCl₂ (0.15 mM). 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 HIF activation by using TransAM™ HIF Kits. The signal intensity for HIF activation in each lot is compared to the signal intensity obtained with extracts from unstimulated COS-7 cells (see figure). After the signals are blanked, the ratio of the signals from stimulated cells over unstimulated cells must be above 4. This ratio may vary depending on the basal level of HIF 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.