

Hep G2 nuclear extract (IL-6 stimulated, 100 ng/ml)

Catalog No: 36107

Species: Human Liver (hepatocellular carcinoma)

Contents: 2 x 100 µg of Hep G2 nuclear extract (IL-6 stimulated, 100 ng/ml) at 2.5 µg/µl

Composition: Nuclear extracts are supplied in lysis buffer consisting 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 b-glycerophosphate, 1 mM DTT and protease inhibitors. The protein content has been determined by a Bradford-based assay.. Cells were cultured in medium supplemented with 100 ng/ml IL-6 (interleukin 6) for 15 minutes at 37°C immediately prior to harvesting.

Quality Control: Each lot has been tested for STAT activation by using TransAM® STAT Family Kits.

Storage and Guarantee: 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. This product is guaranteed for 6 months from date of arrival.