

HeLa whole-cell extract (IL-1 α stimulated)

Catalog No: 40100

Contents: 2 x 125 μ g of HeLa whole-cell extract (IL-1 α stimulated) at 2.50 μ g/ μ l

Composition:

The HeLa whole-cell extracts have been collected in Lysis Buffer after a 30-minute incubation with or without IL-1 α (5 ng/ml). The Lysis Buffer consists of 20 mM Hepes pH 7.5, 350 mM NaCl, 20% glycerol, 1% Igepal-CA630, 1 mM MgCl₂, 0.5 mM EDTA and 0.1 mM EGTA. The protein content has been determined by a Bradford-based assay.

Quality Control:

Each lot has been tested for NF κ B activation by using TransAM™ NF κ B Kits. The signal intensity for NF κ B activation in each lot is compared to the signal intensity obtained with extracts from unstimulated HeLa 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 NF κ B 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.

