

Catalog No: 40050

Species: Human Cervix Adenocarcinoma

Contents: 2 x 100 µg of HeLa nuclear extract at 2.5 µg/µl.

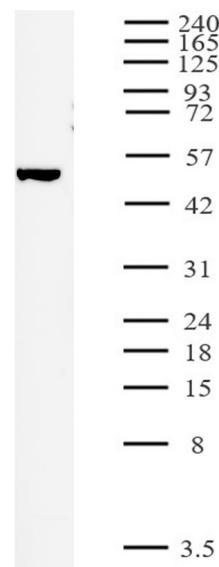
A representative Technical Data Sheet (TDS) is provided here. Please refer to the lot-specific TDS you will receive with your order for the lot-specific buffer contents and protein concentration.

Composition: The HeLa whole-cell extract has been collected in Whole-Cell Lysis Buffer. Whole-Cell Lysis Buffer consists of 25mM Tris-HCl pH 7.5, 150 mM NaCl, 5% glycerol, 1% Igepal-CA630, and 5 mM MgCl₂. The protein content has been determined by a Bradford-based assay.

Quality Control: Extracts have been quality control tested by Western blot.

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.



HeLa whole-cell extract (20 µg/lane)
probed with alpha Tubulin mAb
(1:5,000 dilution).