

HeLa acid extract (Etoposide treated)

Catalog No: 36203

Contents: 100 µg HeLa acid extract (Etoposide treated) at 0.5 mg/ml

Composition:

HeLa acid extracts are supplied in dilution buffer (PBS supplemented with 0.5 mM DTT and 2 mM PMSF). Cells were cultured in medium containing 100 µM etoposide for 6 hours immediately prior to harvesting.

Quality Control:

Acid extracts have been quality control tested by Western blot.

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.