

HeLa acid extract

Catalog No: 36200

Contents: 100 µg of HeLa acid extract at 0.5 µg/µl.

Composition:

Acid extracts are supplied in dilution buffer (PBS supplemented with 0.5 mM DTT and 2 mM PMSF).

Quality Control:

Acid extracts have been quality control tested by Western blot.

Storage:

To ensure stability, extracts should be stored at -80C.

We recommend aliquotting the extracts into single-use fractions and then storing them at -80C. This eliminates repeated freeze/thaw cycles.

Guarantee:

This product is guaranteed for 6 months from date of arrival.