

## HeLa acid extract (Sodium butyrate treated)

---

**Catalog No:** 36202

**Contents:** 100 µg HeLa acid extract (Sodium butyrate 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 5 mM sodium butyrate for 12 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.