

# HeLa Nuclear Extract

**Catalog No.** 17-0001  
**Lot No.** 13059001  
**Pack Size** 250 µg



## Product Description:

HeLa cells are an immortalized human cell line derived from a cervical cancer biopsy. HeLa nuclear extract is an excellent source of transcription factors, histones, chromatin proteins and histone deacetylase (HDAC) activity.

## Formulation:

HeLa nuclear extract at a concentration of 10mg/mL in 20mM HEPES pH 7.5, 400mM NaCl, 1mM EDTA.

## Storage and Stability:

Stable at -80°C for 6 months from date of shipping. For best results, aliquot and avoid repeated freezing and

## Application Notes:

HeLa nuclear extract is useful in a wide variety of experiments, including transcription factor studies, gel shift assays, HDAC assays and as a control for Western blotting.

## References:

This product is for *in vitro* research use only and is not intended for use in humans or animals.