

# HeLa Nuclear Extract



EpiCypher®

**Catalog No.** 17-0001

**Lot No.** 13340002-47

**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 5.33mg/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 freeze/thaw.

## **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:**