

## Ready-to-ChIP HeLa Chromatin

**Catalog No: 53015**

**Format:** 10 rxns

**Quality Control:** 2 vials, each containing 250 µl of HeLa S3 cell sheared chromatin (50 µl @  $1.5 \times 10^6$  cells) for a total of 10 chromatin immunoprecipitation reactions. HeLa S3 cells were grown in suspension, fixed, and the chromatin was prepared using the ChIP-IT® Express protocol for sonication.

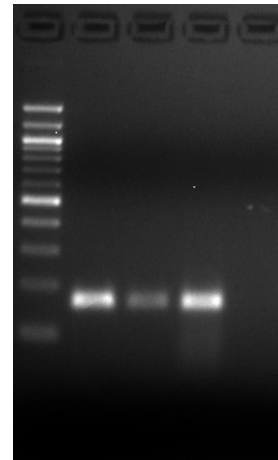
Each lot has been tested for successful chromatin immunoprecipitation with ChIP-IT® Express using a positive control antibody, anti-RNA pol II, and a non-specific IgG as a negative control. The resulting DNA was purified and analyzed by PCR using GAPDH primer control. Positive and negative control antibodies and primers are available in the ChIP-IT Control Kit – Human, Catalog No. 53010.

PCR analysis was performed on DNA isolated through ChIP using the Negative Control IgG (lanes 3 and 7) and the Positive Control antibody (lanes 2 and 6). PCR using the GAPDH primers on DNA immunoprecipitated using the Positive Control antibody generated more product than similar reactions performed on DNA immunoprecipitated using the Negative Control IgG (compare lanes 2 and 3). These results indicate that immunoprecipitation with the Positive Control antibody enriched for GAPDH promoter DNA, and that this enrichment was not the result of non-specific binding of the Positive Control antibody to the chromatin. The input DNA is also assayed by PCR with positive control primers to confirm the presence of the target sequences in the input chromatin (lane 4).

### References:

- Solomon M.J. *et al.* (1988) *Cell* 53(6): 937-47.  
Solomon M.J. and Varshavsky A. (1985) *PNAS USA* 82(19): 6470-4.  
Kuo M.H. and Allis C.D. (1999) *Methods* 19(3): 425-33.  
Weinman A.S. and Farnham P.J. (2002) *Methods* 26: 37-47.  
Caretti G. *et al.* (2003) *J Biological Chem.* 278: 30435-30440.

**Storage and Guarantee:** Store at -80°C. This product is guaranteed stable for 6 months from date of receipt when stored properly.



Chromatin immunoprecipitation using Ready-to-ChIP HeLa Chromatin.

Lane Template Primers  
1 DNA Ladder --  
2 RNA pol II GAPDH  
3 Negative IgG GAPDH  
4 Input DNA GAPDH  
5 H<sub>2</sub>O control GAPDH