

## Ready-to-ChIP NIH/3T3 Chromatin

**Catalog No:** 53021

**Format:** 10 rxns

**Quality Control:** Ready-to-ChIP NIH/3T3 chromatin, consisting of 2 vials each containing 250 µl of NIH/3T3 cell sheared chromatin ( $3 \times 10^4$  cell equivalents per µl,  $1.5 \times 10^7$  cell equivalents total) for a total of 10 chromatin immunoprecipitation reactions. NIH/3T3 cells were grown, fixed, and the chromatin was prepared using the ChIP-IT® Express protocol for sonication.

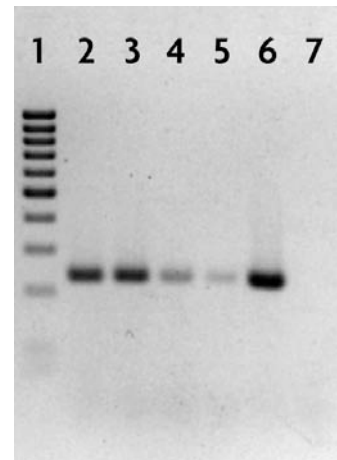
**Quality Control:** Each lot has been tested for successful chromatin immunoprecipitation with ChIP-IT® Express using a positive control antibody, anti-RNA Polymerase II, and a non-specific IgG as a negative control. The resulting DNA was purified and analyzed by PCR using EF1-alpha control primers. Positive and negative control antibodies and primers are available in the ChIP-IT Control Kit - Mouse, Catalog No. 53011.

PCR analysis was performed on DNA isolated from the ChIP reactions using the Positive Control antibody (lanes 2 and 3) and the Negative Control IgG (lanes 4 and 5). PCR using the EF1-α primers on DNA immunoprecipitated using the Positive Control antibody generated reproducibly more product than similar reactions performed on DNA immunoprecipitated using the Negative Control IgG (compare lanes 2 and 3 to lanes 4 and 5). These results indicate that immunoprecipitation with the Positive Control antibody enriched for EF1-α promoter DNA, and that this enrichment was not the result of chromatin binding non-specifically to either the Positive Control antibody or to the beads. The input DNA was also assayed by PCR with positive control primers to confirm the presence of the target sequences in the input chromatin (lane 6).

### 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 NIH/3T3 Chromatin.

- Lane Template Primers  
 1 DNA Ladder --  
 2 RNA pol II EF1-alpha  
 3 RNA pol II EF1-alpha  
 4 Negative IgG EF1-alpha  
 5 Negative IgG EF1-alpha  
 6 Input DNA EF1-alpha  
 7 H<sub>2</sub>O control EF1-alpha