

Ready-to-ChIP K-562 Chromatin

Catalog No: 53020

Format: 10 rxns

Quality Control: Ready-to-ChIP K-562 chromatin, consisting of 2 vials each containing 250 µl of K-562 cell sheared chromatin (3×10^4 cell equivalents per µl, 1.5×10^7 cell equivalents total) for a total of 10 chromatin immunoprecipitation reactions. K-562 cells were grown, fixed, and chromatin was prepared using the ChIP-IT® Express protocol for sonication.

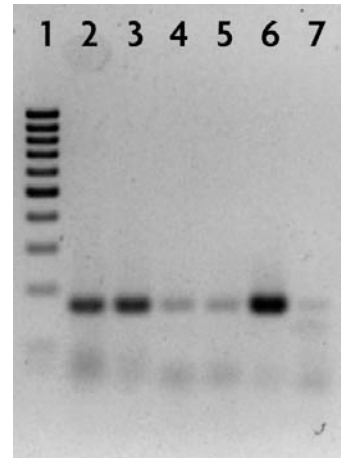
Quality Control: Each lot has been tested for successful chromatin immunoprecipitation with ChIP-IT® Express using a positive control antibody, anti-RNA Polymerase II, and a non-specific IgG as a negative control. The resulting DNA was purified and analyzed by PCR using GAPDH control primers. Positive 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 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 GAPDH 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 GAPDH 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 K-562 Chromatin.

Lane Template Primers
1 DNA Ladder --
2 RNA pol II GAPDH
3 RNA pol II GAPDH
4 Negative IgG GAPDH
5 Negative IgG GAPDH
6 Input DNA GAPDH
7 H₂O control GAPDH