

## Blocker

---

**Catalog No:** 37498

**Format:** 0.1 ml

**Quality Control:** Active Motif's Blocker (part number 103917) is used to reduce non-specific binding during Chromatin Immunoprecipitation (ChIP-Seq) assays. Blocker can be added to the ChIP validated antibody, or mixed with the Protein G beads prior to setting up the immunoprecipitation reaction.

Blocker is provided ready to use at a concentration of 0.5 mg/ml in 100 mM NaCl, 10 mM Tris-HCl and 1 mM EDTA. We suggest using 5  $\mu$ l Blocker with 4  $\mu$ g ChIP-validated antibody in a 30  $\mu$ l reaction volume and incubating for 1 minute at room temperature before adding the antibody/blocker mix to the ChIP reactions. ChIP should then be performed overnight at 4°C with rotation. Alternatively, add 3  $\mu$ l Blocker with 30  $\mu$ l Protein G beads per ChIP and incubate overnight at 4°C with rotation. Then proceed with adding the blocked beads to the immunoprecipitation reaction.

**Storage and Guarantee:** Store at -20°C. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.