

Blocking Reagent AM1

Catalog No: 37496

Format: 0.1 ml

Quality Control: Active Motif's Blocking Reagent AM1 (part number 104891) is used to reduce non-specific binding during Chromatin Immunoprecipitation (ChIP-Seq) assays. Blocking Reagent AM1 can be used to block Protein G beads for use in immunoprecipitation or pre-clearing reactions.

Blocking Reagent AM1 can be used at 3 μ l per 30 μ l Protein G beads. We recommend blocking the Protein G beads overnight at 4°C with rotation.

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.