

ChIP Buffer

Catalog No: 37516

Format: 50 ml

Quality Control: Active Motif's ChIP Buffer (part number 104897) is used in a variety of Active Motif assays to prepare chromatin immunoprecipitation (ChIP) reactions. It is a low SDS (0.1%) buffer to help preserve antibody binding interactions with the target protein during immunoprecipitation.

ChIP Buffer is provided ready to use.

Please note ChIP Buffer is **not** compatible with the ChIP-IT® Express Kits.

Storage and Guarantee: Store at room temperature. 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.