

## 5X IP Low Buffer

---

**Catalog No:** 37511

**Format:** 25 ml

**Quality Control:** The 5X IP Low Buffer (part number 101681) is a component included in Active Motif's Nuclear Complex Co-IP Kit (Catalog No. 54001). The buffer is intended for use in co-immunoprecipitation experiments to study protein-protein complexes. The 5X IP Low buffer is a low stringency buffer that can be supplemented with salt and detergent depending on the complex being studied. Increasing stringency is suitable for stable protein complexes as it reduces background levels. If the strength of the protein complex is not known, we recommend performing experiments evaluating high and low stringency buffers, salt and detergent concentrations to determine the optimal buffer conditions to immunoprecipitate the protein complex.

The 5X IP Low Buffer is provided at a 5X concentration and should be diluted to a 1X concentration using water and/or additional inhibitors, detergents and salt prior to use. For more details, please refer to the product manual for Nuclear Complex Co-IP Kit.

**Storage and Guarantee:** The 5X IP High Buffer is shipped at room temperature and can be stored at 4°C long-term. 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.