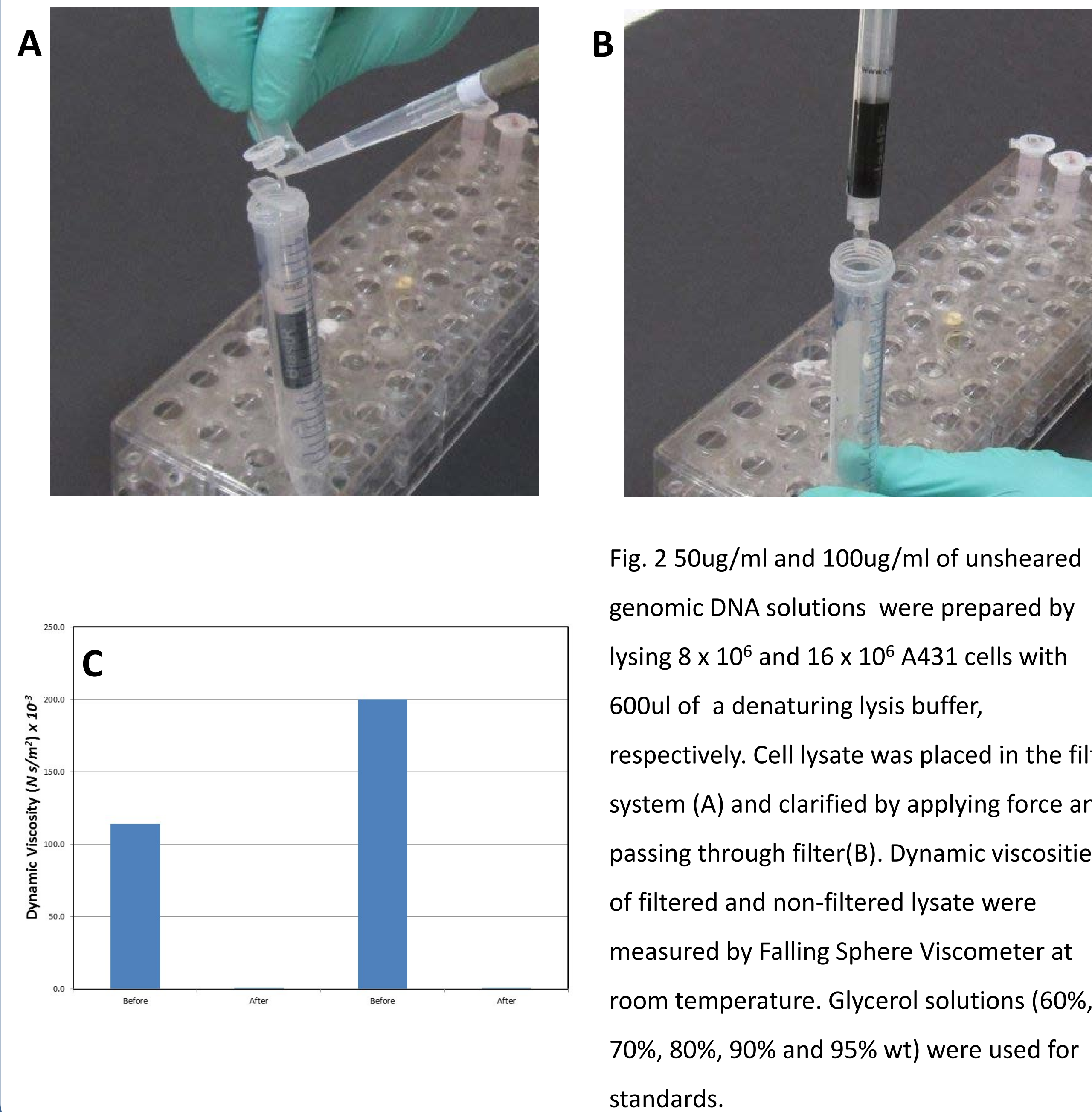


Abstract

The availability of cells in culture provides cell and molecular biologists an invaluable research platform to effectively study signal transduction pathways, protein-expression regulation, and protein-protein interactions. These cells can either be lysed under non-denaturing conditions leading to inefficient recovery of proteins from certain cellular compartments, or with denaturing lysis buffers that efficiently isolates proteins from all cell compartments, but also results in highly viscous genomic DNA (gDNA) contamination. This gDNA contamination can significantly interfere with downstream applications such as western blot (WB) and Immunoprecipitation (IP). However, up until now, no efficient method has been devised to remove a significant portion of this gDNA. Even though the gDNA can be sheared by either passing through a syringe needle or by sonication, small fragments of gDNA can still interfere and/or distort downstream analyses. Additionally, both syringe and sonication methods are highly variable, time consuming, and may even damage protein in the lysate. Moreover, a massive amount of gDNA contamination may obstruct the light transmission in the spectrophotometer, which results in an inaccurate OD reading when obtaining protein concentrations. This study demonstrates a simple and fast method to remove gDNA from the cell lysate completely. A431 cell lysate was purified using this method and more than 98% of gDNA was removed while still recovering 90% of cell lysate without any effect on the protein population. Dynamic viscosities ($N\ s/m^2 \times 10^3$, at room temperature) of the crude cell lysate containing 50ug/ml and 100ug/ml of gDNA were reduced from 114 and 200 to ~ 1 , which is the level of water. Importantly, removal of gDNA resulted in the reliable measurement of the cell lysate protein concentration using a spectrophotometer. We expect that this new method will contribute to the cell biology community by allowing scientist to obtain cell lysates that have a complete protein profile and is free of gDNA contamination. **Contact: tservice@cytoskeleton.com**

2. Filtering reduces the viscosity (gDNA contamination) of cell lysates



Western Blotting (WB)

4. Filtering provides a simple way to obtain high quality cell lysate for WB analysis

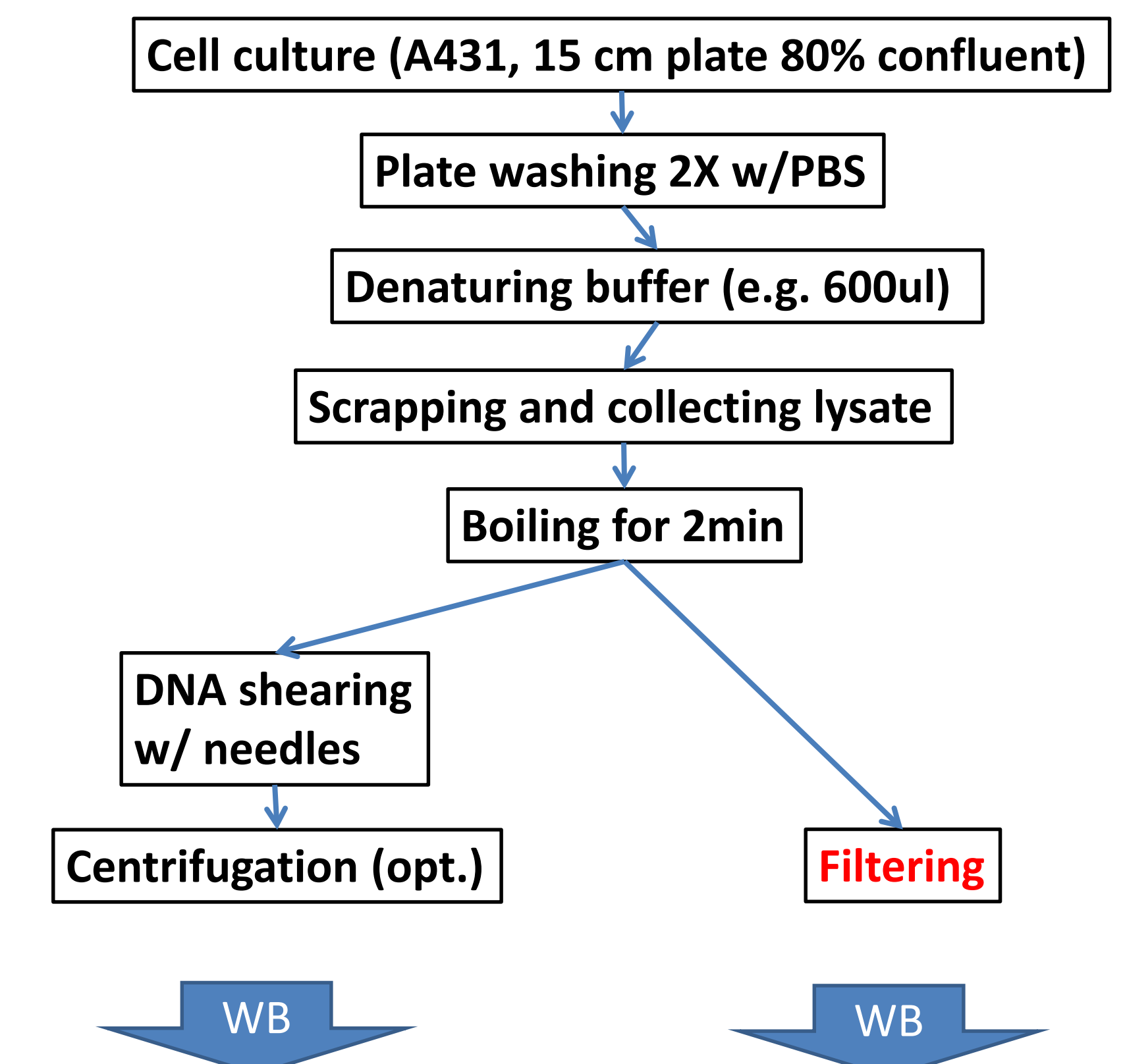


Fig. 4. A comparative schematic diagram between conventional method and new filter method for the preparation of total cell lysate for WB experiment.

Immunoprecipitation (IP)

1. An easy and fast preparation of cell lysate for IP experiments

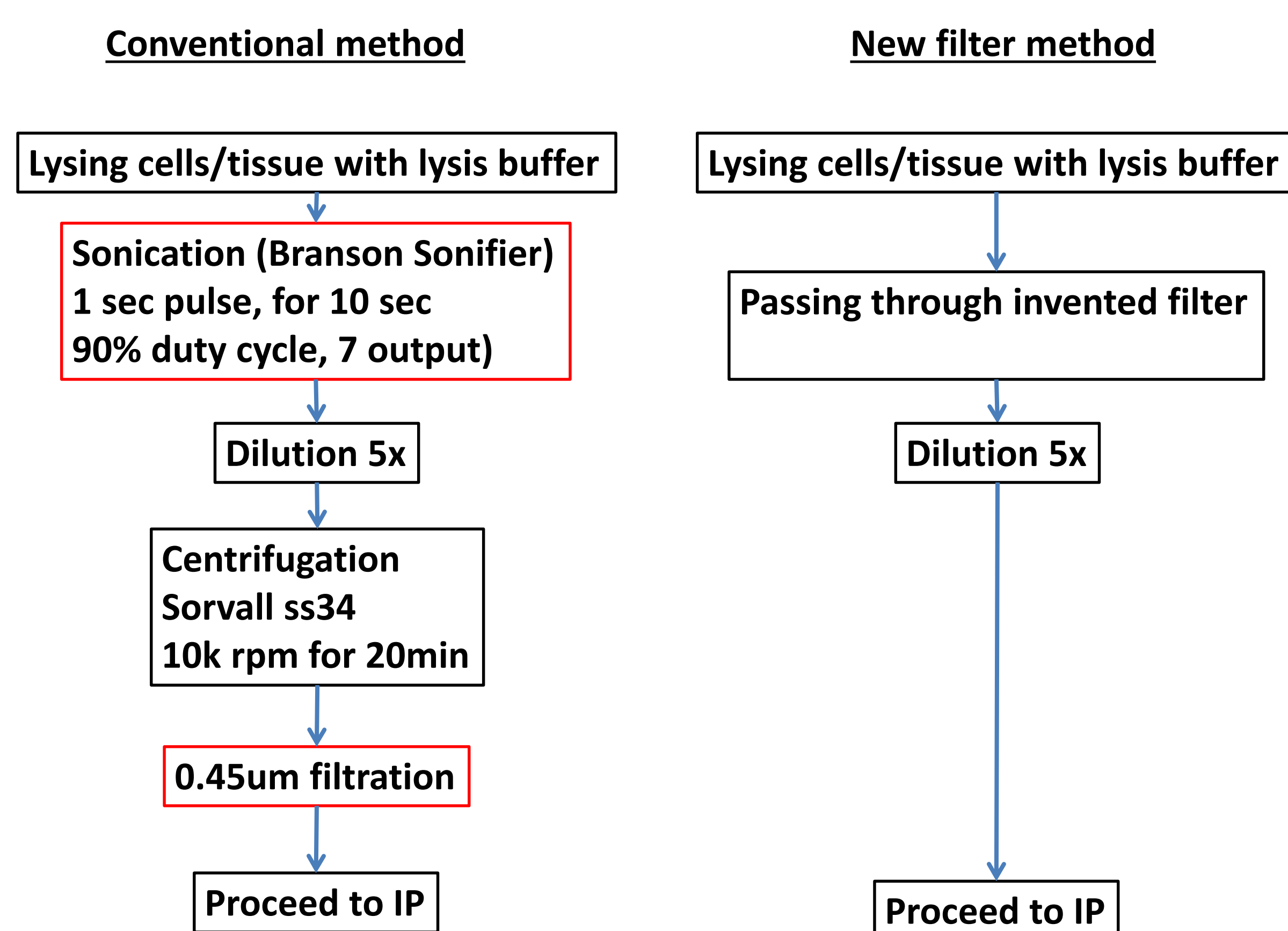
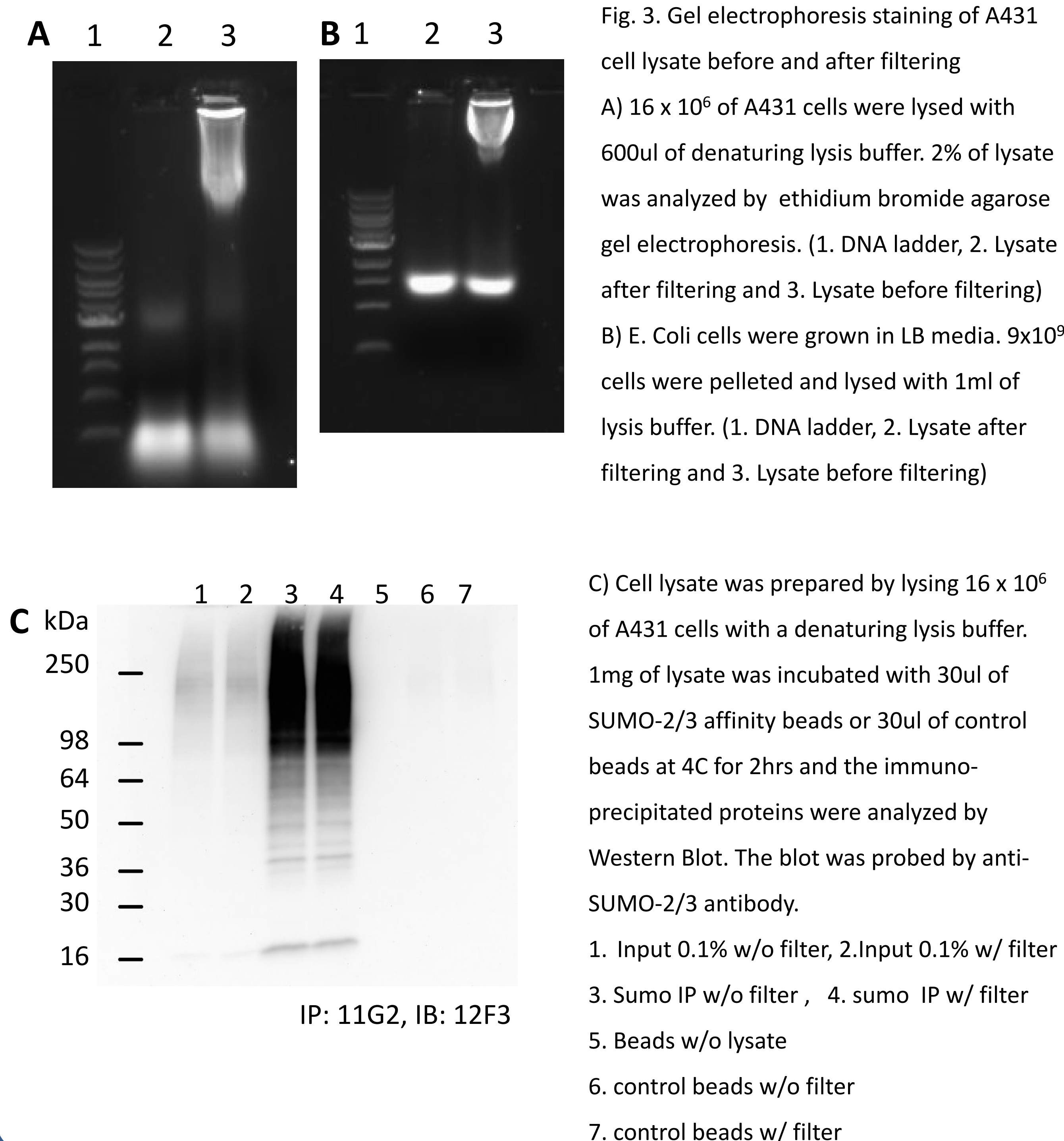
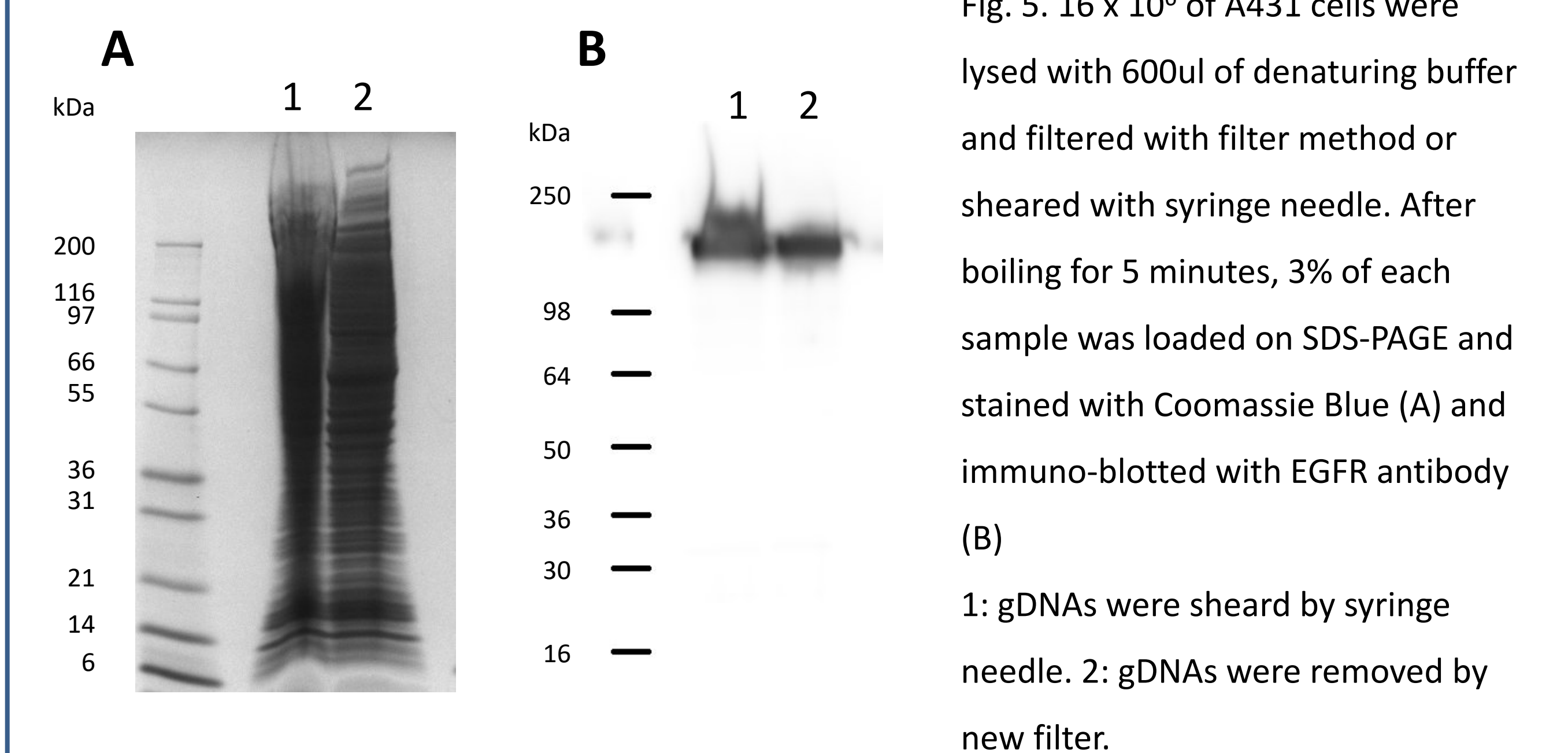


Fig. 1. A comparative schematic diagram between the conventional method and the new filter method for the preparation of total cell lysate for an IP experiment. In order to reduce the gDNA contamination of total cell lysate, the conventional method shears gDNA with either sonication or passing through syringe needle, while the new filter method removes gDNAs with a patented filter system.

3. New filtering method can remove 98% of the gDNA contamination



5. Filtering of gDNA enhances sample quality and analysis by Coomassie staining and western blot



Conclusions

1. New filtering method provides a fast, reliable and convenient method to obtain total cell lysate using denaturing buffers that are convenient for IP and WB experiment
2. This method removes 98% of genomic DNAs from mammalian cell lysate.
3. Dynamic viscosities ($N\ s/m^2 \times 10^3$, at room temperature) of the crude cell lysate containing 50ug/ml and 100ug/ml of gDNAs were reduced from 114 and 200 to ~ 1 , which is the level of water
4. When used with BlastR denaturing buffer, samples can be quantitated using conventional protein quantitation assays