

Mod Spec Histone PTM Quantitation Service

Cells

1. If using cells, prepare 2-5 million cells. For adherent cells, enzymatically detach the cells using trypsin or another enzyme as needed for your specific cell type. If cells grow in suspension, just transfer cells in growth media to a conical tube for pelleting.
2. Spin at 800 x *g* for 5 minutes to pellet cells and then remove the supernatant.
3. Wash cells once in PBS and transfer to a 1.7 ml microfuge tube.
4. Spin at 800 x *g* for 5 minutes to pellet cells.
5. Remove supernatant with a pipette and freeze the cell pellet on dry ice. Store at -80°C.

Tissues

1. If using tissues, isolate 25 to 100 mg of tissue.
2. Place the tissue in a 1.7 ml microfuge tube and freeze by placing the tube in liquid nitrogen or by placing the tube on dry ice. Store at -80°C.