

## 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.

### 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.
- Ship cells on dry ice. Fill out Active Motif's **Sample Submission Form**, include a hard copy with your samples and ship according to the instructions on the Sample Submission Form.

**North America** 877 222 9543

**Europe** +32 (0)2 653 0001

**Japan** +81 (0)3 5225 3638

**www.activemotif.com**

**North America** 877 222 9543

**Europe** +32 (0)2 653 0001

**Japan** +81 (0)3 5225 3638

**www.activemotif.com**

**Mod Spec Sample Prep RevA\_092020**