

Sample Preparation Protocol

When sending your samples for our analytical pipeline please send cell-free supernatant. Should you not know how to achieve that please prepare them according to this protocol.

1. Centrifuge an appropriate volume of fresh cell culture sample for 5-10 min at 200 x g.
2. Filter supernatant with a 0.2 µm filter into an appropriate tube/container.
3. Store your samples at your appropriate temperature, e.g. -20 °C and ship accordingly e.g. on dry ice.

Important information:

- You should prepare your samples according to this sample preparation protocol before sending them to us.
- Your samples have to be free of bacterial and viral contamination.
- Your samples have to be non-hazardous and classified as Biosafety Level 1.
- Please attach the sample declaration sheet to your samples.
- If different storage conditions are necessary or desired, please contact us
- For accurate information on how and where to send your samples please contact us directly