

Please ship your samples to:

Sartorius Xell GmbH
FAO Dr. Tim Steffens
Waldweg 21
33758 Schloss Holte-Stukenbrock
Germany
Opening hours: 8.00am - 5.00pm

Date

Dear Sartorius Xell GmbH,

we hereby declare our samples to be (check appropriate boxes):

- free of cells and debris
 free of bacterial and viral contamination

The Biosafety Level of our samples is: _____

- non-hazardous
 free of toxic | radioactive material

or:

- our samples contain (e.g., antibiotics): _____
 prepared according to Xell's sample preparation protocol

or:

- prepared differently (specify): _____

Kind regards,

Company; Name; Signature

Important information:

- Please attach this declaration to your samples and/or send a scan in advance
- Please include the Sartorius order number with your samples. You will find this on the order confirmation
- You should prepare your samples according to our sample preparation protocol
- Samples should be sent on dry ice/wet ice (as appropriate)!
- Any further questions, please contact a member of our sales team