



## Chitozen Q&A

### Q/ How do I recognize the side of the coverslip that is chitosan-coated?

The chitosan-coated side does not touch the microscope slide. It allows to avoid frictions that would damage the coating. Thus, it is the external side of the coverslip that is the one you must use, as indicated on the following scheme. Stickers will also help you in finding the coated side.



### Q/ Am I obliged to centrifuge the slides?

You are not obliged to centrifuge the slides.

Idylle guarantees a good adhesion of the following bacteria

- with centrifugation: Escherichia coli
- without centrifugation: Bacillus subtilis, Pseudomonas fluorescens

**Q/ Does the culture medium affect bacteria adherence?**

No. It rather has an effect on the growth speed.

**Q/ Do you have any recommendations about the flux?**

Tâm Mignot and colleagues apply a flux of 1,5 mL/min, which is fine when bacteria adhere correctly to the chitosane.

We recommend not to apply a flux higher than 5 mL/min.

**Q/ Which bacteria have been validated with Chitozen?**

Escherichia coli, Vibrio cholerae, Myxococcus xanthus, Mycobacterium smegmatis, Bacillus subtilis, Pseudomonas aeruginosa and Pseudomonas fluorescens.

September 2, 2022.