



SAMPLING OF *P. JIROVECI* IN THE SURROUNDING AIR OF PATIENTS WITH PNEUMOCYSTIS PNEUMONIA

Saint Louis Hospital, APHP, France

/ CONTEXT

Pneumocystis jirovecii is responsible for a disease affecting immunodepressed patients: *Pneumocystis pneumonia* (PCP). Airborne transmission of *Pneumocystis* has been demonstrated in animal models and is highly probable in humans. However, information concerning burdens of *Pneumocystis jirovecii* in exhaled air from infected patients is lacking due to the technical limitations of the current sampling methods. This fungus is indeed not cultivable, so impaction method based on the cultivability of microorganisms cannot be used. Also, the difficulties of extraction from a solid support or the lengthy conditions of sampling (filtration technique) led us to consider alternative methods that would be suitable for both sampling and quantifying *P. jirovecii* in air. Our objective is to evaluate *P. jirovecii* air diffusion in patients with *Pneumocystis pneumonia* with Coriolis air sampler sampling combined with RT-PCR assays [1].

/ PROTOCOL

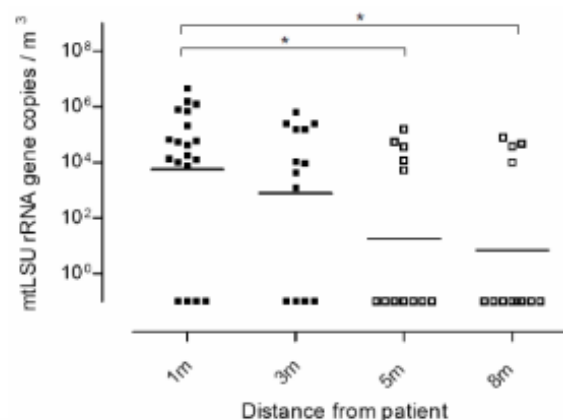
- Collection of air samples in each patient room with the Coriolis® μ air sampler.
- Collection of 1 air sample of 1.5 m³ (5 min) at a distance of 1 m from patient's head + air samples inside and outside the room: at the door entrance and in the corridor (3m, 5m and 8m from the patient).
- Centrifugation of the liquid sample at 1300 g for 10 min; removal of the supernatant to leave a 1 mL pellet; storage at +4°C.
- Extraction of DNA from 200 μ L of sample pellets using the QIAamp DNA Mini Kit (Qiagen).
- Performance of a real-time quantitative PCR assay targeting the mitochondrial large subunit (mtLSU) rRNA gene of *P. jirovecii* (Applied Biosystems 7500 PCR system).

/ RESULTS

33/56 positive air samples : the more the distance from the patient is low, the more the *P.jirovecii* DNA is detected at:

1m - 15/19 samples (78.9%), 3m - 9/13 (69.2%), 5m - 5/12 (41.7%), 8m - 4/12 (33.3%)

This study provides the first quantitative data on the spread of *P. jirovecii* in exhaled air from infected patients.



Number of *P. Jirovecii* mtLSUrRNA gene copies per m³ of air sample inside (■) and outside (□) the room. *P<0.05; Dunn's post test.v

/ CONCLUSION

The Coriolis® samples large volumes of air during a short period and concentrates the airborne microorganisms into a small volume of liquid. This was crucial for efficient capture and detection of *Pneumocystis*. This facilitates epidemiological investigations with non-cultivable microorganisms (like *Pneumocystis* sp.) in conjunction with real-time PCR amplification of DNA.

This study leads the way for initiating a quantitative risk assessment for airborne transmission of *P. jirovecii* and other airborne infectious agents.

Authorised distributor



In Australia:
For customer service, call 1300-735-292
To fax an order, use 1800-067-639
To email an order, ordersau@thermofisher.com

In New Zealand:
For customer service, call 0800-933-966
To fax an order, use 0800-329-246
To email an order, ordersnz@thermofisher.com



CNIM group