

Fluorescence analysis of reactive oxygen species (ROS) generated by six isolates of *Aspergillus fumigatus*

U-C. Hipler¹, U. Wollina¹, D. Denning²

¹Department of Dermatology, Friedrich-Schiller-Universität, Jena, Germany ²Department of Microbiology, University of Manchester, UK

- Detection of intracellular H₂O₂
- LOD of 5 nM H₂O₂ in 200 µL
- LOQ of > 10⁷ cells per mL

Introduction

Reactive oxygen species (ROS) are essential intermediates in oxidative metabolism. Nonetheless, when generated in excess, ROS can damage cells by peroxidizing lipids and disrupting structural proteins, enzymes and nucleic acids. Excess ROS are generated during a variety of cell stresses, including ischemia/reperfusion, exposure to ionizing and ultraviolet radiation and/or inflammation. ROS may contribute to inflammation and tissue damage.

The processes leading to ROS generation can be monitored using luminescence analysis or fluorescence methods. Intracellular ROS generation in cells can be investigated using 2',7'-dichlorofluoresceindiacetate (DCFH-DA), which is an established compound to detect and quantify intracellularly produced H₂O₂.

The conversion of non-fluorescent DCFH-DA to the highly fluorescent compound, 2',7'-dichlorofluorescein (DCF), happens in several steps. First, DCFH-DA is transported across the cell membrane and deacetylated by esterases to form the non-fluorescent 2',7'-dichlorohydrofluorescein (DCFH). This compound is trapped inside of the cells. Next, DCFH is converted to DCF through the action of peroxide, which is generated by the presence of peroxidase.

Aspergillus species are of interest in the pathogenesis of several dermatological diseases. It is uncertain whether *Aspergillus* itself may generate ROS and therefore actively induce tissue damage. The present study investigates whether *Aspergillus* species are capable of producing ROS by themselves and if there are differences between the several strains.

Materials & Methods

Six isolates of *Aspergillus fumigatus* (AF 65, AF 71, AF 72, AF 91, AF 210, AF 294) cultured 5 weeks on Sabouraud-Glucose-Agar (BAG) were investigated.

After addition of isotonic NaCl solution and centrifugation with 100 rpm for 10 min, the blastospore concentration could be estimated by counting in CASY 1 (Schärfesystem).

These cell suspensions with concentrations of 10⁵ to 10⁷ cells/mL were measured on a filter-based BMG LABTECH microplate reader using 100 µL of fungal cell suspension after incubation with 100 µL DCFHDA (0.4 nM) for each single test. To eliminate LBS induced effects, polymyxin B (3 mg/mL) was added to all experimental suspensions.

Each measurement was done at least sixteen times in duplicate for calculation of the mean and the standard error of the mean.

Results & Discussion

The ability of various *Aspergillus* species to generate ROS was investigated.

For all fungal cells, a linear increasing fluorescence activity could be observed depending on the incubation time with DCFH-DA. By using a calibration curve, the measured fluorescence signals were converted to H₂O₂ concentrations and one example is given for AF 71 with different incubation periods (Fig. 1). Because of the small sized *Aspergillus fumigatus* (2-3 µm), detectable fluorescence was observed only at concentrations >10⁵ cells/mL.

The ROS generation showed a linear and direct proportional dependence on cell numbers and the results were reproducible on 3 different days with a fixed incubation period of 2.5 hours at 37°C (Fig. 2). The highest value could be found at concentrations of 10⁷ cells/mL.

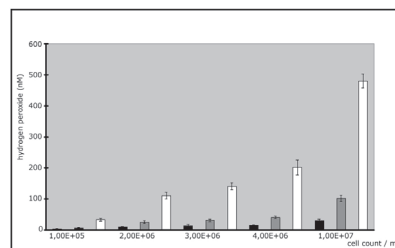


Fig. 1: H₂O₂ production of AF 71 depends on incubation period and cell number.

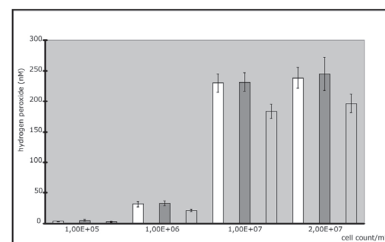


Fig. 2: H₂O₂ production of AF 71 depends on cell number with 4 hr incubation time at 37°C on three separate days.



The isolates of *A. fumigatus* AF 91 and AF 72 are resistant against itraconazole (antifungal agent) and AF 65 is resistant to amphotericin B (antifungal agent). We investigated whether or not there are connections between the resistance and the ROS generation. Interisolate differences could be found (Fig. 3).

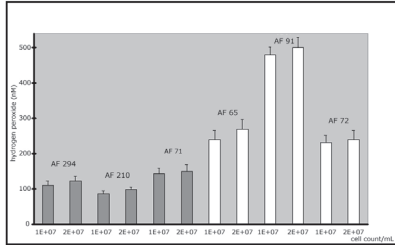


Fig. 3: H₂O₂ production of several strains of *Aspergillus fumigatus* with 3 hr incubation time at 37°C. The last three strains are resistant against antifungal agents.

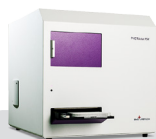
Conclusion

The morphological event of fungi, usually acknowledged as a major factor of virulence, is associated with increased intracellular ROS formation, which are most likely secreted. Together with phospholipases, ROS are capable of destroying the host cell membranes. This process may contribute to the invasiveness of *A. fumigatus* and to the inflammatory response of the host.

Clearly, the pathogenicity of *Aspergillus* species is a function of a multitude of parameters working together in a sequential and cooperative manner to establish infection.

It has been shown by several authors, that water as well as superoxide, hydrogen peroxide and OH-radicals can be generated in the course of the mitochondrial electron transport process. Fungi also possess the normal and the alternative pathway of electron transport. The interesting connection between the monovalent oxygen reduction and the energy conservation in isolated chloroplasts was described. The hydrogen peroxide formation at the phosphorylation points I and II is disposed by a cytochrome c oxidation process.

In this study, the method of ROS fluorescence measurement was utilized on different unstimulated *Aspergillus* species for the first time. There were linear correlations found between ROS levels and blastospore concentrations. A pathophysiological meaning of the released oxygen metabolites as an additional factor of virulence in the complicated system of inflammatory reactions, which was also estimated in *Saccharomyces cerevisiae*, is hence not to be excluded.



PHERAstar® FSX

*The PHERAstar FSX is the newest PHERAstar reader.



CLARIOstar®



Omega Series