

# FRET sensor assesses ATP levels in living plants experiencing low oxygen conditions

Stephan Wagner<sup>1,2</sup>, Philippe Fuchs<sup>1,2</sup>, Thomas Nietzel<sup>1,2</sup>, Marlene Elsässer<sup>1,2</sup>, and Markus Schwarzländer<sup>1,2</sup>

<sup>1</sup>Institute of Plant Biology and Biotechnology, University of Münster, Germany <sup>2</sup>Institute of Crop Science and Resource Conservation (INRES), University of Bonn, Germany

- The genetically encoded sensor ATeam monitors (sub)cellular MgATP levels *in vivo* and in real-time
- The CLARIOstar® microplate reader can record MgATP dynamics in the plant cytosol, e.g. at oxygen shortage
- Low oxygen stress results in depletion of cellular MgATP pools in living plants

## Introduction

To survive as sessile organisms, plants need to constantly adapt their metabolism to their environment. Flooding of plants is widespread and has severe metabolic consequences as it limits the cellular supply with oxygen to drive respiration, and thus production of ATP. Submergence acclimation typically involves drastic alterations in metabolism to circumvent anoxia and to maintain primary metabolism and ATP supply. We have recently established the use of a genetically encoded protein sensor [ATeam1.03-nD/nA, see Ref. 1] to assess MgATP dynamics in the model plant *Arabidopsis thaliana*. This system allows for reliable analyses of changes in MgATP *in vivo* and in real-time and provides a new handle to investigate the energetic consequences of low oxygen for the plant cell.

## Assay Principle

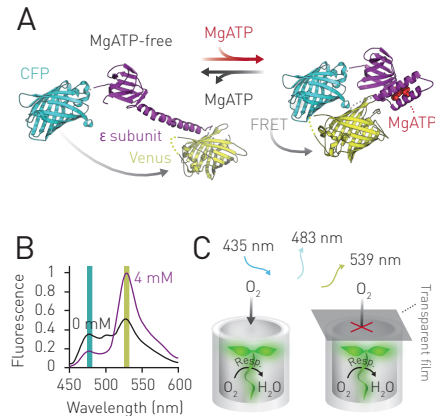
ATeam is a genetically encoded Förster resonance energy transfer (FRET)-based protein sensor that reversibly binds MgATP<sup>1</sup>. MgATP binding induces a structural change of the MgATP-binding domain and results in more efficient FRET from the blue fluorophore mseCFP (CFP) to the yellow fluorophore cp173-mVenus (Venus). (Figure 1A). High Venus/CFP fluorescence emission ratios thus indicate high MgATP levels and vice versa (Figure 1B). ATeam can be expressed stably, constitutively and at high levels in *Arabidopsis* plants and low oxygen stress can be imposed by submerging plants while simultaneously restricting their oxygen supply (Figure 1C).

## Materials & Methods

- CLARIOstar microplate reader (BMG LABTECH)
- Transparent Microtest Plate 96 Well,F (Sarstedt #82.1581)
- Transparent film for qPCR (VWR #391-1258)
- Assay medium (10 mM MES, pH 5.8, 10 mM CaCl<sub>2</sub>, 10 mM MgCl<sub>2</sub>, 5 mM KCl)

### Experimental procedure

*Arabidopsis* seedlings were submerged in assay medium on a transparent 96 well microtest plate. Wells were either left unsealed (control condition allowing for resupply of the medium with atmospheric oxygen) or sealed with a transparent film to restrict resupply of the assay medium with oxygen. The ATeam fluorescence was recorded while - except from the excitation flashes - the plate was kept in the dark to suppress oxygen evolution through photosynthesis.



**Fig. 1:** **(A)** Structural model of ATeam. The MgATP-binding domain [ε subunit of ATP synthase] is flanked by CFP and Venus. The structural transition from MgATP-free to MgATP-bound form increases the FRET efficiency between the two fluorophores. **(B)** Binding of MgATP at increasing concentrations between 0 and 4 mM leads to a fluorescence emission decrease of the FRET donor CFP (at 483 nm) and a fluorescence emission increase of the FRET acceptor Venus (at 539 nm). **(C)** *Arabidopsis* seedlings stably expressing ATeam in the cytosol were exposed to low oxygen stress by restricting oxygen supply to plants submerged in assay medium. Residual oxygen in sealed wells is gradually consumed through active respiration (Resp.). Submerged seedlings in unsealed wells served as control. Figure adapted from Ref. 2.

### Instrument settings

Optic settings	Fluorescence intensity, plate mode kinetic, top optic
	Number of multichromatics: 2
Monochromator settings	CFP [gain 1250] Excitation 435-20 nm Dichroic 459.5 Emission 483-18
	Venus [gain 1200] Excitation 435-20 nm Dichroic 488.8 Emission 539-13
Well scan	Orbital averaging with a 3 mm diameter
General settings	Number of flashes per cycle: 50
	Settling time: 0.2 s
Kinetic settings	Number of cycles: 800
	Cycle time: 180 s
Shaking	10 s double orbital shaking at 400 rpm before each cycle



## Results & Discussion

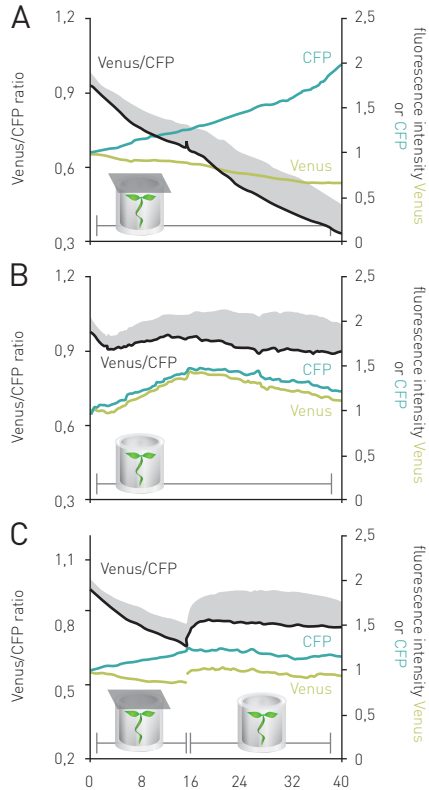
Arabidopsis seedlings exposed to low oxygen stress showed gradually decreasing Venus/CFP ratios in the cytosol (Figure 2A), while the values remained almost constant in seedlings that were submerged in assay medium but not sealed to allow for oxygen re-supply, except for an initial decrease immediately following seedling immersion (Figure 2B). This suggests that MgATP supply to the cytosol under oxygen-limiting conditions is not able to balance cytosolic ATP consumption. Unsealing of wells after 16 h led to a recovery of MgATP levels through re-oxygenation of seedlings, demonstrating the reversibility of ATeam readings (Figure 2C). Individual emission intensity plots of Venus and CFP highlight a key advantage of ratiometric FRET-based protein sensors over intensimetric sensors and/or dyes: while in Figure 2B individual fluorophore intensities vary [e.g., through sample movement], the Venus/CFP ratio remains relatively stable. A true difference in MgATP binding as seen in Figure 2A is, in contrast, characterized through the antiparallel behaviour of both fluorophore emission intensities as a result of *bona fide* FRET.

## Conclusion

The CLARIOstar microplate reader allows to read the fluorescence of genetically encoded protein sensors (e.g., ATeam1.03-nD/nA) expressed in plants exposed to naturally occurring stress conditions [e.g., hypoxia as a hallmark of submergence stress].

## References

1. Imamura, H et al. (2009) Visualization of ATP levels inside single living cells with fluorescence resonance energy transfer-based genetically encoded indicators. *Proc. Natl. Acad. Sci. USA* **106**: 15651-15656.
2. De Col, V et al. (2017) ATP sensing in living plant cells reveals tissue gradients and stress dynamics of energy physiology. *eLife* **6**: e26770.



**Fig. 2:** Plant cytosolic MgATP dynamics under low oxygen stress. Ten-day-old Arabidopsis seedlings expressing cytosolic ATeam and grown *in vitro* were submerged in assay medium. Three seedlings in each well were excited at  $435\pm 10$  nm and the emission of CFP ( $483\pm 9$  nm) and Venus ( $539\pm 6.5$  nm) was read using well multichromatics. Wells in **(A)** were sealed with a transparent film to limit oxygen resupply; wells in **(B)** were left unsealed as a control and wells in **(C)** were first sealed and then un-sealed after 16 h to allow re-oxygenation of seedlings. Each panel shows the normalized Venus/CFP ratio (black, primary axis) averaged from 12-16 wells. Error bars (grey) are SD. Additionally, each panel shows the normalized emission intensities of Venus (yellow) and CFP (cyan) averaged from the same wells (secondary axis). Figure adapted from Ref. 2.



PHERAstar® FSX



CLARIOstar®



Omega Series