

Determination of Maximal Chlorophyll Fluorescence Using A Multiphase Single Flash of Sub-Saturating Intensity

S. D. Loriaux¹, R. A. Burns¹, J. M. Welles¹, D. K. McDermitt¹, and B. Genty²

¹LI-COR Inc., 4421 Superior St., P. O. Box 4425, Lincoln, NE, USA

²CEA/Cadarache DEVM-Laboratoire d'Écophysiologie Moléculaire des Plantes, UMR 6191-CNRS-CEA-Univ. Aix Marseille II, 13108 Saint Paul lez Durance, Cedex, France.

INTRODUCTION

- It is necessary to determine maximal fluorescence (F_m') in order to estimate the effective quantum efficiency of Photosystem II (Φ_{PSII}) and the rate of electron transport (ETR) from chlorophyll fluorescence measurements (Genty *et al.*, 1989).
- F_m' is commonly measured using a single saturating, multi-turnover flash of 400-1200 ms duration to reduce the Q_A -PQ pool. In many conditions, full reduction of the Q_A -PQ pool cannot be achieved using such a protocol, especially in field plants.
- Here we present a 'multiphase' single flash method (MSF) to estimate the true maximal fluorescence level in these conditions.

MATERIALS & METHODS

- With light emitting diode (LED) technology, "saturating" light can be supplied in several different ways, which are described in Figure 1.
- LI-6400 Portable Photosynthesis Systems with 6400-40 Leaf Chamber Fluorometers (LI-COR, Lincoln, NE, USA) were used to compare the three methods. Chamber conditions were maintained at ambient levels of CO_2 , relative humidity, and temperature.
- Φ_{PSII} was calculated as: $1 - F_s/F_m'$, where F_s is steady state fluorescence and F_m' corresponds to maximal fluorescence. ETR was calculated as: $\Phi_{PSII} \cdot P \cdot Q \cdot \alpha_{Lef}$. The fraction of absorbed quanta used by PSII, f , was assumed to be 0.5; incident photon flux density, Q , was measured with an internal PAR sensor; and maize leaf absorbance, α , was measured as 0.85.

THE MULTIPLE FLASH TRAIN SOLUTION

- Markgraf & Berry, 1990 and Earl & Ennahli, 2004, proposed an approach using a multiple flash train (FT) of variable light intensities (Q) to estimate the true F_m' (Figure 1).
- Table 1 shows the rectangular single flash (RSF) method under-estimates F_m' compared to the FT method at all light intensities in field-grown maize leaves.
- However, the FT method requires multiple fluorescence measurements at different light pulse intensities, with wait times between each measurement, to obtain one valid estimate of F_m' , Φ_{PSII} and ETR, which takes about 10 minutes. This is time-consuming and laborious if many measurements are to be made, and is especially difficult under rapidly changing environmental and physiological conditions.

THE NEW "MULTIPHASE" SINGLE FLASH SOLUTION

- The MSF method uses a ramp of rapidly varying light intensity to generate F_m' values over a range of light intensities (B, Figure 1).
- Total duration of one MSF measurement is only about one second.

Validation with DBMIB

- As a first step to evaluate the MSF method, we vacuum infiltrated leaves of *P. sativum* with 150 μM DBMIB to restrict turnover of the PQ pool and provide conditions in which nearly full saturation could be assured.

- The fluorescence response during the ramp phase (Panel II Inset, Figure 2) has a hyperbolic shape. The y-intercept of a plot of the fluorescence response as a function of $1/Q$ (Panel II, Figure 2) is used to estimate the maximal fluorescence at infinite Q .
- F_m' measured by the RSF method was within 1% of the extrapolated F_m' value measured by the MSF method (Panel II, Figure 2). Thus it should be possible to use the MSF method to derive F_m' in conditions where saturation by the traditional RSF method cannot be achieved.

Validation in field-grown *Z. mays*

- Figure 3 illustrates the new MSF method in field-grown maize. In this case, the traditional RSF method gave a 17% lower estimate of F_m' than the extrapolated value predicted by the MSF method.
- Table 2 and Figure 4 show that the FT and MSF methods gave similar estimates of F_m' , Φ_{PSII} and ETR at all light levels.
- Figure 5 shows that ETR calculated by both the new MSF and FT methods are proportional to gross CO_2 assimilation at all irradiances, as expected for C_4 plant metabolism; however, ETR estimated using the traditional RSF method increasingly underestimates results from gas exchange as irradiance increases.
- In maize, under-estimation of F_m' by the traditional RSF method can lead to miscalculation of Φ_{PSII} and ETR by as much as -41% at high light intensities. The new multiphase single flash (MSF) method resolves this problem.

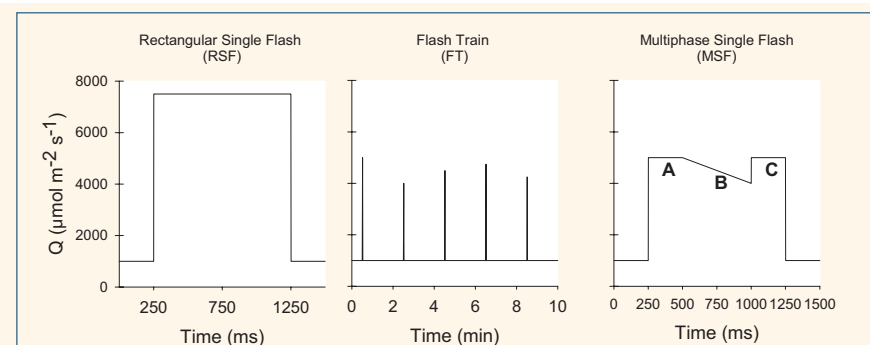


Figure 1. Methods to determine maximal fluorescence (F_m').

Rectangular Single Flash method (RSF): a saturating multi-turnover flash (Q) of 400-1200 ms duration. **Flash Train method (FT):** applies five rectangular flashes of various Q in random order, separated by two minutes. **Multiphase Single Flash method (MSF):** (A) high, nearly saturating Q for 250 ms to reduce Q_A -PQ pool; (B) ramp of declining Q for 500 ms; (C) return to the initial high Q for 250 ms to check for flash-induced non-photochemical quenching (qN). F_m' values from phase B of the MSF method and each flash of the FT method are regressed against $10^4/Q$ and extrapolated to estimate the maximal fluorescence at infinite flash intensity.

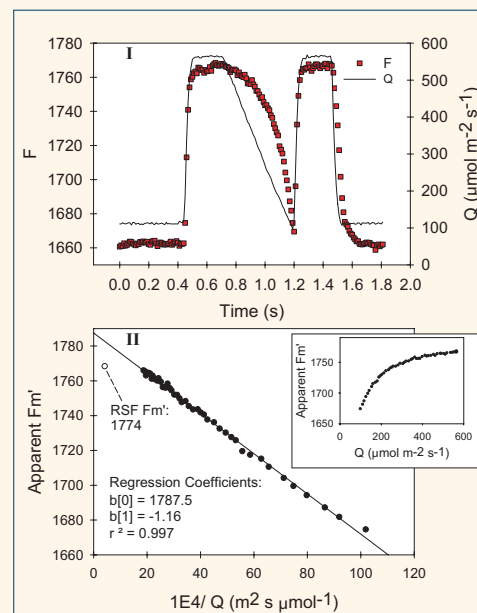


Figure 2. Panel I. Fluorescence response to the MSF protocol in a typical leaf of greenhouse-grown *P. sativum* vacuum infiltrated with DBMIB.

Panel II. Linear regression of F_m' versus $10^4/Q$ obtained during phase B, Panel I.

Inset: Fluorescence response to Q during the ramp.

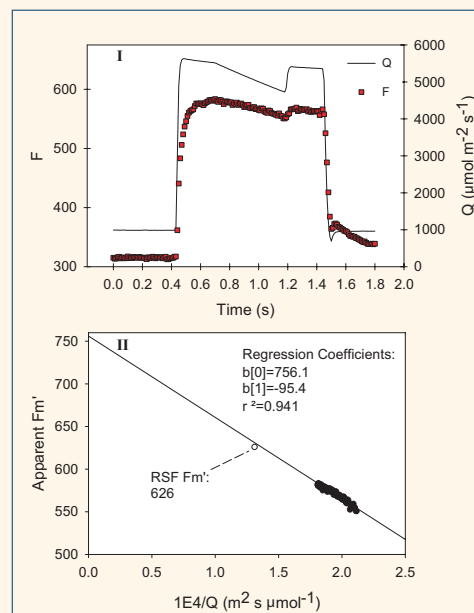


Figure 3. Panel I. Typical fluorescence response to the MSF protocol in a field-grown maize leaf at an incident photon flux density of $1,000 \mu mol m^{-2} s^{-1}$. Panel II. Linear regression of F_m' versus $10^4/Q$ using the ramp data.

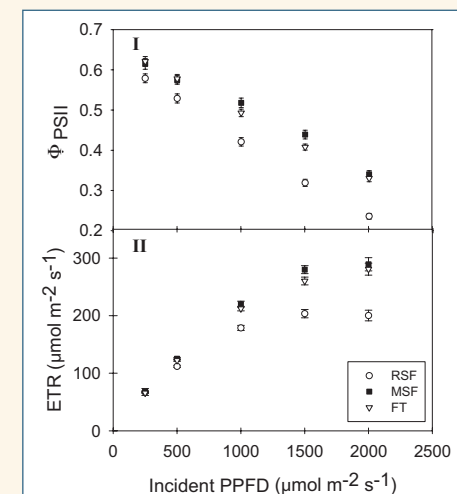


Figure 4. Panel I. Φ_{PSII} measured in field-grown maize at growth stage V-5 as a function of light intensity. Panel II. PSII electron transport rate (ETR) computed from the data in Panel I versus light intensity. Each data point is the mean of 9 to 14 observations. Error bars indicate standard errors of the means.

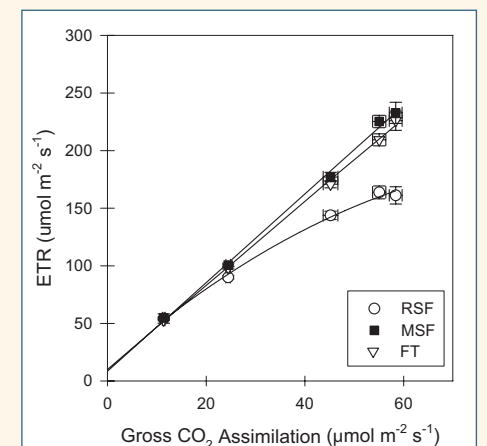


Figure 5. Comparison of ETR calculated using the RSF, FT and MSF methods to gross CO_2 assimilation rate measured from gas exchange in field-grown maize at growth stage V-5.

Table 1. Comparison of F_m' , Φ_{PSII} , and ETR in field-grown maize leaves measured by RSF and FT methods. Data are computed as $[(RSF - FT)/FT] \times 100\%$ for each of the variables on each of n leaves.

Incident PPFD ($\mu mol m^{-2} s^{-1}$)	n	F_m'	Φ_{PSII} , ETR
250	9	-19**	-8**
500	11	-16**	-10**
1000	14	-20**	-19**
1500	13	-22**	-28**
2000	14	-16**	-41**

** Significant at $\alpha 0.01$

Table 2. Comparison of F_m' , Φ_{PSII} , and ETR in field-grown maize leaves measured by FT and MSF methods. Data are computed as $[(MSF - FT)/FT] \times 100\%$ for each of the variables on each of n leaves.

Incident PPFD ($\mu mol m^{-2} s^{-1}$)	n	F_m'	Φ_{PSII} , ETR
250	9	2 ^{ns}	3 ^{ns}
500	11	6 ^{ns}	2 ^{ns}
1000	14	4 ^{ns}	4 ^{ns}
1500	13	4*	8*
2000	14	6 ^{ns}	4 ^{ns}

^{ns} Significant at $\alpha 0.05$

*Significant at $\alpha 0.05$

REFERENCES

- Genty, B., J.-M. Briantais, and N. R. Baker. 1989. The relationship between the quantum yield of photosynthetic electron transport and quenching of chlorophyll fluorescence. *Biochim. Biophys. Acta* 990: 87-92.
- Earl, H. J. and S. Ennahli. 2004. Estimating photosynthetic electron transport via chlorophyll fluorometry without photosystem II light saturation. *Photosynth. Res.* 82:177-186.
- Markgraf, T. and J. Berry. 1990. Measurement of photochemical and non-photochemical quenching: correction for turnover of PS2 during steady-state photosynthesis. *In: M. Baltscheffsky (ed.), Curr. Res. Photosynth.* IV:279-282.

CONCLUSIONS

- The new multiphase single flash method can be used to derive true estimates of F_m' , Φ_{PSII} and ETR in a flash time duration of approximately 1s.
- The traditional rectangular saturating flash method leads to large under-estimates of F_m' , Φ_{PSII} and ETR, especially at higher light intensities.
- For the first time, this new approach allows both rapid and valid determination of the rate of photosynthetic electron transport in conditions where F_m' is difficult to saturate, or when leaf physiological and environmental status is changing rapidly (e.g. field conditions).

©2006. LI-COR is a registered trademark of LI-COR, Inc. All other trademarks belong to their respective owners. LI-COR is an ISO 9001 registered company.

This work was supported by LI-COR Biosciences, www.licor.com

LI-COR
Biosciences