

Fluorometric determination of extracellular enzyme activities in peat using a BMG LABTECH microplate reader

Bonnett, S.A.F., Leah, R., and Maltby, E.

Institute for Sustainable Water, Integrated Management and Ecosystem Research, University of Liverpool, Merseyside, UK

- Fluorescence method to measure extracellular enzyme activities in peat
- Methylumbelliferone (MUF) and derivatives used as substrates
- Fast measurement in 96-well format using a BMG LABTECH microplate reader

Introduction

Peatland ecosystems represent a significant global store of carbon (C) due to the historic accumulation of organic matter resulting from suppressed microbial de-composition. Decomposition is limited by extracellular enzymes that are produced by the plant and microbial communities and function independently of the microbial community. The storage of carbon in peatland ecosystems is therefore regulated in part by the environmental factors that determine extracellular enzyme activities. Understanding the environmental controls and processes that regulate extracellular enzyme activities is therefore critical in determining the fate of carbon in peatland ecosystems.

Here we present the results from a method developed to measure potential extracellular enzyme activities in peat using a BMG LABTECH microplate reader. The assay is based on the use of fluorogenic methylumbelliferyl (MUF) substrates that fluoresce upon enzymatic cleavage allowing the amount of product to be measured.

Materials & Methods

- Black 96-well microplates, F-bottom, Greiner Bio-One
- MUF and MUF substrates, Sigma-Aldrich
- BMG LABTECH microplate reader

Peat samples were collected from Langdon moor, Durham, UK in order to determine the concentration of methylumbelliferyl (MUF) substrates to be added to peat solutions to achieve enzyme 'active-site' saturation and the optimal assay incubation length.

The hydrolytic enzymes β -glucosidase, cellobiohydrolase, and N-acetylglucosaminidase (or chitinase) were determined using MUF artificial substrates 3 [see Table 1]. For each peat replicate sample, ten cm³ of peat was placed in a 50 mL centrifuge vial and deionised water added up to the 50 mL mark. The vial was shaken by hand for 30 seconds and mixed using a vortex for a further 30 seconds. Using a sawn-off pipette tip, 0.75 mL of peat slurry was placed into a 1.7 mL centrifuge vial. To the vial, 0.75 mL of MUF substrate was added and samples were incubated at field temperature for one hour. Samples were centrifuged for 5 minutes at 12,000 rpm and 300 μ L of supernatant transferred to a well in a plate. Fluorescence was determined on a BMG LABTECH microplate reader at 330 nm excitation and 450 nm emission wavelength. Enzyme activities were determined from the fluorescence units using a standard calibration curve of methylumbelliferone (MUF) and expressed as rates of MUF production (μ mol MUF per g⁻¹ dry peat weight per min⁻¹). Fluorescence quenching is a potentially interfering process which decreases the

intensity of the fluorescence emission and occurs in water containing peat-derived compounds. The standard calibration curve accounted for quenching by dissolving the MUF standard in 150 μ L of centrifuged peat slurry for each peat replicate.

Table 1: Enzymes and enzyme MUF substrates.

Enzyme	MUF substrate	Reaction
Cellobiohydrolase	MUF- β -D-cellobioside	Cellulose polymers to dimers
β -glucosidase	MUF- β -D-glucopyranoside	Cellulose dimers to monomers
Chitinase	MUF-N-acetyl- β -D-glucosaminide	Chitin to acetylglucosamine

The assay above was performed with substrate concentrations varying from 0 to 500 μ M in order to determine the optimal substrate concentration. Using the optimal substrate concentrations, samples were then incubated for varying lengths of time to determine the optimal assay incubation length.

Results & Discussion

Figure 1 shows the effect of increasing substrate concentration on the activity of cellobiohydrolase. Activity followed Michaelis-Menten kinetics and reached saturation at approximately 200 μ M of substrate.

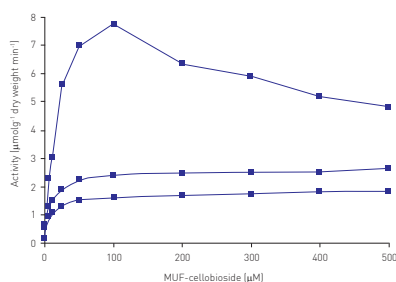


Fig. 1: Effect of substrate concentration on cellobiohydrolase activity (n = 3).

One of the replicates exhibited substrate inhibition at 100 μ M of substrate resulting in a relative standard deviation of 85 % at 100 μ M of substrate. This is mainly due to spatial variation of enzyme activity within the block of peat being analysed, NOT analytical variation.

Figure 2 shows the effect of increasing substrate concentration on the activity of β -glucosidase. Activity followed

Michaelis-Menten kinetics and reached saturation at approximately 200 μM of substrate. One of the replicates exhibited substrate inhibition at 100 μM of substrate resulting in a relative standard deviation of 91 % at 100 μM of substrate.

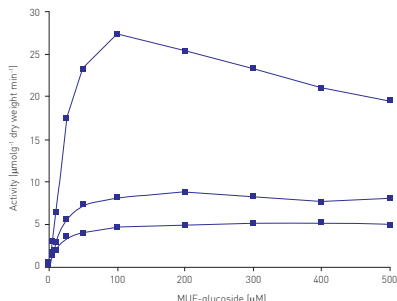


Fig. 2: Effect of substrate concentration on β -glucosidase activity ($n = 3$).

Figure 3 shows the effect of increasing substrate concentration on the activity of chitinase. Activity followed Michaelis-Menten kinetics and reached saturation at approximately 300 μM of substrate.

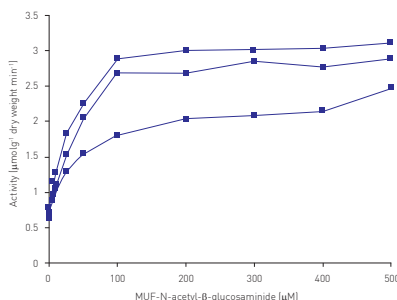


Fig. 3: Effect of substrate concentration on chitinase activity ($n = 3$).

Figure 4 shows the significant quenching effect of peat on fluorescence.

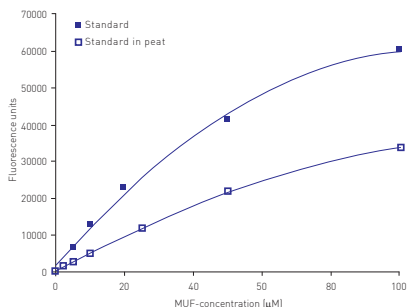


Fig. 4: The quenching effect of peat on fluorescence.

Time course incubations performed using optimal substrate concentrations for cellobiohydrolase, β -glucosidase and chitinase showed that one hour was the maximum amount of time that reaction rates remained linear (data not shown).

Conclusion

The extracellular enzymes β -glucosidase, cellobio-hydrolase and N-acetylglucosaminidase reached substrate saturation at 200 μM , 200 μM and 300 μM of MUF respectively. Measured enzyme activities were linear for up to 120 minutes although 60 minutes duration of incubation was chosen for future applications to minimise variation between replicates. One replicate exhibited substrate inhibition for the β -glucosidase and cellobio-hydrolase assays.

The results show that potential extracellular enzyme activities can be determined in peat at low cost and within a short period of time using a BMG LABTECH microplate reader.

