

# Luminescent ABEL® Antioxidant Test Kit with PHOLASIN® for Vitamin C Type Antioxidants



Bernd Hipler, BMG LABTECH GmbH, Germany  
Jan Knight, Knight Scientific Limited, UK



## 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 stresses on cells including ischemia / reperfusion, exposure to ionizing and ultraviolet radiation and/or inflammation. Reactive oxygen species may contribute to inflammation and tissue damage. As an example, consider the fate of the all important polyunsaturated lipids in a cell membrane. Attack by a free radical on one of the carbon to carbon double bonds can lead to the formation of the corresponding lipid peroxy radical and the initiation of a chain reaction in which the final products, lipid peroxides, no longer maintain the integrity of the cell wall.

This process, lipid peroxidation, may be initiated by the hydroxyl radical OH<sup>•</sup> in a reaction in which iron, as

Fe<sup>2+</sup>, is involved. In biological systems much attention is focused on the two free radicals, superoxide anion and hydroxyl radical.

The generation processes of ROS can be monitored using luminescence. Because of the very weak native luminescence of ROS, both luminol and lucigenin have been used in the past to give measurable signals. However, the luminescence of these two substances is very small compared to that of Pholasin®, which allows precise analysis on very small samples.

Pholasin® is the photoprotein of the marine rock boring bioluminescent mollusc, Pholas dactylus and the Common Piddock. Pholasin® does not glow on its own, but can be switched on by free radicals and other reactive oxygen species. The capacity of samples to scavenge free radicals and other oxidants can be monitored by means of the described test kit that contains Pholasin®.

## Results and Discussions

A white 96-well plate loaded with the assay components and different samples was placed in the LUMIstar and the reagent injector was primed with solution B. The assay was done as described above. If Pholasin® is present when the superoxide is generated, light will be emitted.

If there are antioxidants in the sample capable of scavenging superoxide, then these antioxidants will compete with Pholasin® for the superoxide and less light will be detected. Controls containing no sample are run with each assay.

The luminescence measurements with the ABEL® Antioxidant Test using different substances with unknown antioxidant capacity are shown in Figure 3.

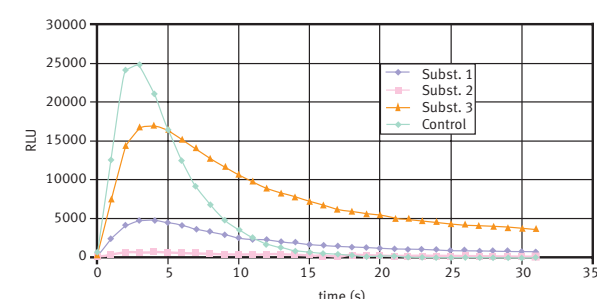


Fig. 3: Measurement of 3 different substances with ABEL® Antioxidant Test: Superoxide Assay

These substances were prepared from natural plant extracts. Compared with the control (100%) scavenging of superoxide could be detected. The inhibition amounted 32%, 80% and 97%, respectively. It is possible to quantify the antioxidant capacity of unknown substances by means of the ascorbate standard curve (Fig. 4).

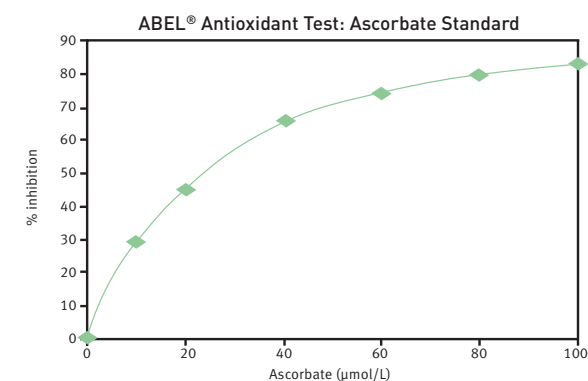


Fig. 4: Ascorbate Equivalent Antioxidant Units

The inhibition of 80% corresponds to about 77 μmol/L in the well. The Ascorbate Equivalent Antioxidant Units in the sample (10 μL) is obtained by multiplying the x value (i.e. 77 μmol/L) by the dilution factor (20) to obtain 1.5 mmol/L.

## Materials and Methods

All materials were obtained through normal distribution channels from the manufacturer stated.

- LUMIstar, BMG LABTECH, Offenburg, Germany



Fig. 1: The LUMIstar microplate reader

- Test kit KSL-ABEL-21 M2, Knight Scientific Limited (15 Wolseley Close Business Park, Plymouth, PL2 3BY, UK) e-mail: Knightscientific@aol.com, Web site: www.knightscientific.com
- 96-well plate, white, BMG LABTECH, Offenburg, Germany
- Samples of unknown antioxidant capacity

Although the LUMIstar was utilized for this application note, BMG LABTECH's FLUOstar, POLARstar or NOVostar could also have been used for luminescence measurements. In addition, consumables such as pipette tips and tubes were used as needed from various manufacturers.

### Protocol for Antioxidant Test for Superoxide

Superoxide is generated instantaneously when Solution B is added to Solution A. By means of the reagent injector Solution B is injected into the microplate containing Solution A together with Pholasin®, with and without a 10 μL sample of fluid of unknown antioxidant capacity.

Into a microplate well add the following :

- 25 μL Assay Buffer (for controls) or 15 μL Assay Buffer + 10 μL sample
- 50 μL Antioxidant Pholasin®
- 100 μL Solution A

The assay is initiated when 25 μL of Solution B is injected into the microplate well, which must be in the light measuring position.

The assay is carried out using the following parameters:

- Gain = 200, well mode
- No. of intervals = 32
- Meas. Interval Time = 1 s
- Total Meas. Time = 32 s
- Start injection = 0 s

### Standard curve for Ascorbate

Pipette into a microplate well :

- 50 μL Pholasin
- 5 μL Reconstitution and Assay buffer (R&A)
- 100 μL Solution A

Using the 2 built-in reagent injectors of the LUMIstar the dilution series of ascorbate can be generated automatically. Injector 1 is programmed with the volumes of 1 mmol/L ascorbate solution and injector 2 with the volumes of R&A buffer as shown in Table 1.

Table 1: Volumes of ascorbate and R&A buffer added to microplate for recording a standard curve

1 mmol/L ascorbate (μL)	R&A Buffer (μL)	Ascorbate (μmol/L) in 200 μL
0	20	0
2	18	10
4	16	20
8	12	40
12	8	60
16	4	80
20	0	100

The corresponding ascorbate standard curve is shown below (Fig. 2).

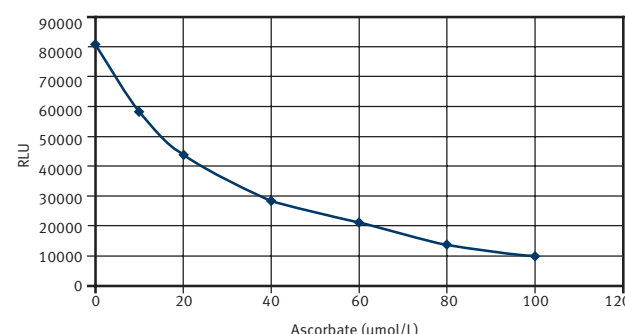


Fig. 2: Ascorbate standard curve

## Conclusion

The ABEL® Antioxidant test kit with Pholasin® can be used successfully for measurements of substances with unknown antioxidant capacity. Using the luminometer LUMIstar the test can be done very fast and with a high precision and reproducibility.

Germany: BMG LABTECH GmbH Tel: +49 781 96968-0  
Australia: BMG LABTECH Pty. Ltd. Tel: +61 3 59734744  
China: BMG LABTECH Co. Ltd. Tel: +86 10 6424063  
France: BMG LABTECH SARL Tel: +33 1 48862020  
UK: BMG LABTECH Ltd. Tel: +44 1296 336650  
USA: BMG LABTECH Inc. Tel: +1 919 806 8526  
www.bmglabtech.com info@bmglabtech.com