

ORAC Assay Performed on the POLARstar Omega and PHERAstar FS microplate reader

Franka Ganske and Marjan Orban, BMG LABTECH, Offenburg, Germany

Application Note 197

Rev. 07/2009



- Antioxidants are able to neutralize Reactive Oxygen Species
- Antioxidant capacity of fruit juices determined with the ORAC assay in 96 and 384-well format
- MARS data analysis software with predefined templates for quick ORAC evaluation

Introduction

In all oxygen consuming cells, metabolism and oxidative stress generate several intermediates and byproducts that are collectively known as reactive oxygen species (ROS). ROS are necessary intermediates in the human body, but they are also involved in the aging process and in the development of many degenerative diseases, including cancer, heart disease, Alzheimer's and Parkinson's. ROS are dangerous to cellular structures and functional molecules (i.e. DNA, proteins, lipids) as they act as strong oxidizing agents or free radicals. Biological antioxidants are able to dispose of ROS; however, they are not completely effective in eliminating all of the free radicals, oxygen ions and peroxides that can do damage to the body. Furthermore, ROS can be generated from exposure to other external sources such as cigarette smoke, pollutants, chemicals and environmental toxins.

One standardized method for determining the antioxidant capacity of a substance is the ORAC (oxygen radical absorbance capacity) assay.¹ The ORAC assay is based upon the inhibition of the peroxy-radical-induced oxidation initiated by thermal decomposition of azo-compounds such as [2,2'-azobis(2-methylpropionamide) dihydrochloride (AAPH)].²

Herein we describe the application of the ORAC assay on a POLARstar Omega and PHERAstar FS (Fig. 1) using Trolox® (a water-soluble analogue of vitamin E) as a standard substance. Different fruit juices were measured for their ORAC value.

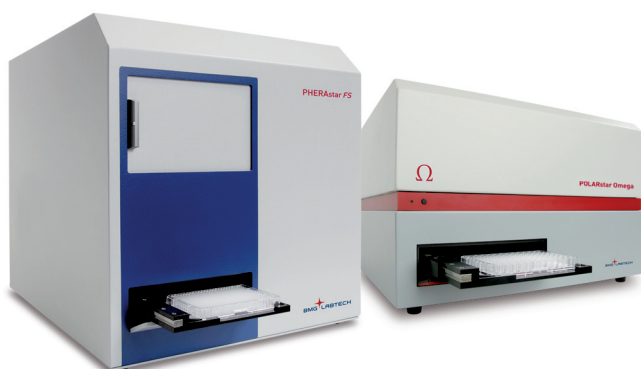


Fig. 1: BMG LABTECH's multidetection microplate readers PHERAstar FS and POLARstar Omega

Assay Principle

Over time ROS, generated from the thermal decomposition of AAPH, will quench the signal from the fluorescent probe fluorescein. The subsequent addition of an antioxidant produces a more stable fluorescence signal, with signal stability depending on the antioxidant's capacity. Data points are summarized over the time by the MARS data analysis software. This is then compared to the standard, Trolox®, and is expressed as micromoles of Trolox® equivalents (TE per gram or per milliliter of sample (μmole of TE/g or μmole of TE/mL). Please refer to Application Note 148 to see the assay principle graphically.³

Materials and Methods

- Black 96-well and 384-well plates from Greiner (#655076 and #781076)
- Fluorescein sodium, 6-hydroxy-2,5,7,8-tetra-methylchroman-2-carboxylic acid (Trolox®), L (+)-ascorbic acid, Epicatechin gallate, [2,2'-azobis(2-methylpropionamide) dihydrochloride (AAPH)] were obtained from Sigma-Aldrich
- Plate sealer, BMG LABTECH, Aylesbury, UK, Cat. No. 77400-05
- POLARstar Omega, BMG LABTECH, Offenburg, Germany
- PHERAstar FS, BMG LABTECH, Offenburg, Germany

Eight different fruit juices were obtained from a local supermarket. All juices were indicated to contain 100 % fruit juice except for the peach and pear juice which were labelled to contain 50 % fruit juice.

Test Protocol

Different dilutions of Trolox® (12.5 μM to 200 μM) and fruit juices were prepared in phosphate buffer (10 mM, pH 7.4). All solutions should be prepared fresh daily.

In every working well of a 96-well plate the following was pipetted in triplicate. For a 384-well plate half of the volumes were used:

- 150 μL of a 10 nM Fluorescein solution (all wells)
- +25 μL Trolox® dilution (standards)
- +25 μL sample dilution (samples)
- +25 μL phosphate buffer (blank)

The microplates were sealed followed by an incubation for 30 min at 37°C in a THERMOstar microplate incubator. Alternatively, incubation can be performed in the microplate reader.

After incubation, fluorescence measurements (ex. 485 nm, em. 520 nm) were taken every 90 sec to determine the background signal. After 3 cycles, 25 μL of 240 mM AAPH was injected using the onboard injectors. Alternatively, AAPH can also be added manually with a multi-channel-pipette. This has to be done as quickly as possible since the ROS-generator displays immediate activity after addition. The test was resumed and fluorescence intensity measure-

ments were taken up to 120 minutes. An instrument setting overview is given in application note 148.³

Results and Discussion

Fruit juices have ingredients that can quench the fluorescence signal. This usually depends on the dilution used (Fig. 2).

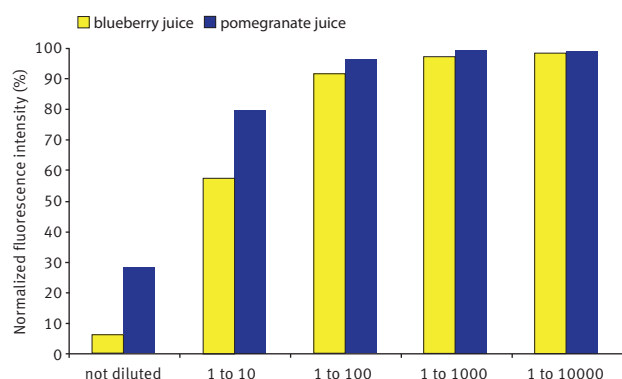


Fig. 2: Normalized fluorescence values of 10 nM fluorescein (prepared in phosphate buffered saline) in presence of 15 % blueberry and pomegranate juice (for different dilutions).

Figure 2 shows that not diluted blueberry juice quenches the fluorescence signal below 10 % of the expected value whereas diluting leads to nearly no quenching. The degree of quenching varies between the different fruit juices and can cause problems in ORAC evaluation when the data is not normalized.

Signal curves taken from the MARS data analysis software in figure 3 show the fluorescence decrease of fluorescein after addition of the ROS generator AAPH over time. The area under the signal curve is depending on the concentration of the antioxidant standard substance (Fig. 3).

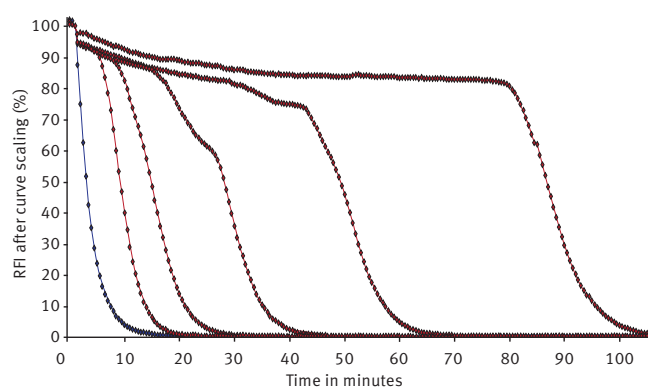


Fig. 3: Signal curves for different Trolox® concentrations (red graphs) and a blank without Trolox (blue graph) recorded on the PHERAstar FS in 384-well format. The curves were normalized to 100 %. The 100 % value is the maximum value that is obtained directly after injection of AAPH.

The area under the curve of samples is needed to calculate the Trolox® equivalents. In case of quenching ingredients, that are still active even in high dilutions, all signal curves should be normalized. This can be easily done with the MARS data analysis software using the “Curve Scaling” feature. Applying Curve Scaling will allow for optimal area under the curve calculation.

The Trolox® standard curve allows for back calculation of samples. In Table 1 a summary of all tested fruit juices is given.

Table 1: ORAC values in µmol TE/mL not diluted fruit juice. Assay was measured on the POLARstar Omega and PHERAstar FS in 96-well and 384-well format. The data shows an average of three independent measurements.

	POLARstar Omega		PHERAstar FS	
	96-well	384-well	96-well	384-well
Blueberry juice	66.9	66.0	56.5	62.5
Pomegranate juice	34.1	31.6	26.0	31.3
Sallow thorn juice	64.7	75.0	69.5	67.9
Apple juice	6.5	5.0	6.2	5.5
Grape juice	12.2	11.9	14.6	11.8
Orange juice	16.3	15.5	15.5	15.3
Peach juice	8.8	7.3	7.4	7.5
Pear juice	9.1	7.5	6.8	6.2

From the tested fruit juices the sawallow thorn and blueberry ones show the highest ORAC values. Please notice that the values may differ when compared to similar measurements as juice ingredients are different from manufacturer to manufacturer.

Conclusion

It was shown that the ORAC assay can be easily performed on the POLARstar Omega and the PHERAstar FS using the onboard injectors. The results obtained in both plate formats, 96-well and 384-well, are similar. Further downscaling of the total volume may be possible in combination with the PHERAstar FS.

MARS data analysis software comes with every Omega and PHERAstar reader and allows automated evaluation. Predefined templates in combination with curve scaling features give ORAC results quickly and easily.

References

1. Cao, G., Alessio, H. M., Cutler, R. G. (1993) Oxygen-radical absorbance capacity assay for antioxidants. *Free Radical Biol. Med.* **14**, 303-311.
2. Glazer, A. N. (1990) Phycoerythrin Fluorescence-Based Assay for Reactive Oxygen Species. *Methods Enzymol.* **186**, 161-168.
3. BMG LABTECH Application note 148 (2006): <http://www.bmglabtech.com/application-notes/fluorescence-intensity/orac-148.cfm>

Germany: BMG LABTECH GmbH **Tel: +49 781 96968-0**

Australia: BMG LABTECH Pty. Ltd. **Tel: +61 3 59734744**
France: BMG LABTECH SARL **Tel: +33 1 48 86 20 20**
Japan: BMG LABTECH JAPAN Ltd. **Tel: +81 48 647 7217**
UK: BMG LABTECH Ltd. **Tel: +44 1296 336650**
USA: BMG LABTECH Inc. **Tel: +1 919 806 1735**

Internet: www.bmglabtech.com **info@bmglabtech.com**