

# DPPH Antioxidant Assay Kit

## Technical Manual

Technical Manual (Japanese version) is available at <http://www.dojindo.co.jp/manual/d678.pdf>

### General Information

Recent findings suggest that a decline in internal antioxidant capacity causes the onset of various diseases and health impairment. Consequently, interest in antioxidant rich foods has been recently increasing. Shimamura et al. improved the 2,2-diphenyl-1-picrylhydrazyl (DPPH) assay, a procedure for evaluating antioxidant capacities, and reported it may be applicable as a standard method to evaluate the antioxidant capacity of antioxidants.<sup>1)</sup>

The DPPH Antioxidant Assay Kit is based on the DPPH assay improved by Shimamura and enables quick and easy measurements of the antioxidant capacity of a sample. Using this kit, the antioxidant capacity is expressed as the Trolox equivalent antioxidant capacity (TEAC), a value calculated from the IC<sub>50</sub> of the antioxidant and the IC<sub>50</sub> of Trolox. The DPPH Reagent and Trolox Standard included in this kit only need to be dissolved prior to use. Moreover, the protocol for this kit supports a microplate assay format for the simultaneous analysis of multiple samples

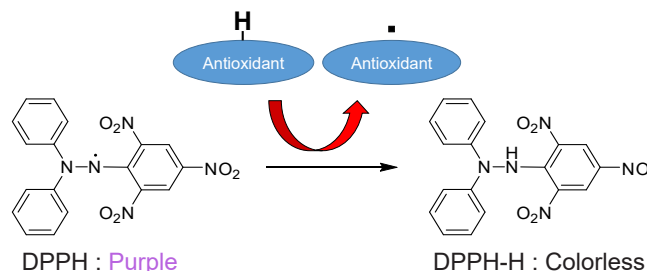


Figure 1. Principle of the DPPH Antioxidant Assay Kit

### Kit Contents

	100 tests	500 tests
DPPH Reagent	× 1	× 5
Trolox Standard	1 mg × 1	1 mg × 5
Assay Buffer	11 mL × 1	55 mL × 1

### Storage Conditions

- Store at 0–5 °C

### Required Equipment and Materials

- Ethanol (99.5%)
- 10 mL measuring flask
- Microplate reader (equipped with a filter around 517 nm)
- 96-well Microplate
- Incubator (25°C)
- Micropipettes
- 20–200 µL Multichannel pipette
- Ultrasonic cleaner

### Precautions

- Equilibrate reagents to room temperature prior to use.
- Each tube of DPPH Reagent and Trolox Standard is sufficient for 100 tests.
- Please note that the contents may be on the tube wall or on the ceiling of the cap.
- Analyzing samples and standards in triplicate is recommended for accuracy.

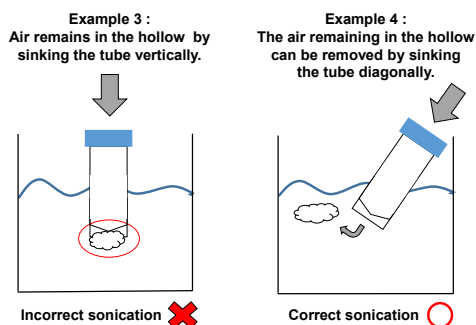
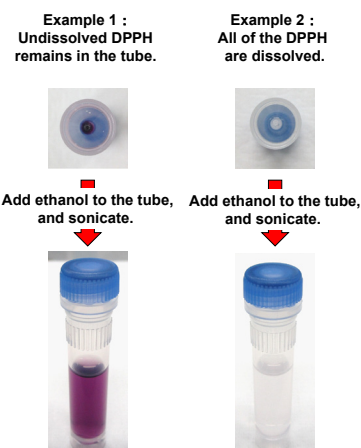
### Preparation of Solutions

#### Preparation of the DPPH working solution

1. Add approximately 1 mL of ethanol to a tube of DPPH Reagent and sonicate for 60 seconds.
2. Transfer all of the solution prepared in step 1 to a 10 mL measuring flask.
3. Add another aliquot of approx. 1 mL of ethanol to the tube from step 1 and sonicate for 60 seconds.

4. Check the color of the solution and if the solution has a purple color as shown in the images on the right, repeat steps 2 and 3 until no color is observed.

- \* Please use sonication to dissolve the DPPH Reagent because it is difficult to dissolve.
- \* Undissolved DPPH may result in the variation of data. Please make sure that all of the DPPH is dissolved.
- \* If air remains in the hollow of the tube, it is hard to dissolve the DPPH Reagent completely by sonication. Please remove the air by sinking the tube of the DPPH Reagent diagonally into a ultrasonic cleaner.



5. Make up to a final volume of 10 mL with ethanol.
- \* Prepare the DPPH working solution fresh each day.

## Preliminary Experiment

### Preparation of the Trolox Standard solution

1. Add approximately 1 mL of ethanol to the Trolox Standard tube and completely dissolve the contents by vortexing or sonication.
2. Transfer all of the solution prepared in step 1 to a 10 mL measuring flask and add ethanol to 10 mL.
3. Dilute the 100 µg/mL Trolox Standard solution prepared in step 2 with ethanol to make 80, 60, 40, and 0 µg/mL solutions (Figure 2).

\* Prepare the Trolox Standard solution fresh each day.

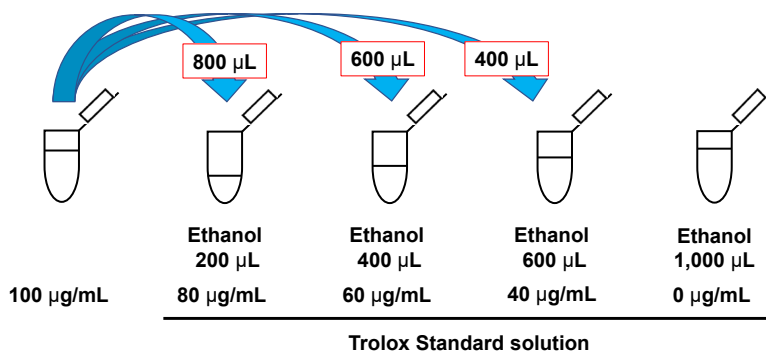


Figure 2. Preparation of the Trolox Standard solutions

- TEAC is calculated based on the IC<sub>50</sub> values for the Trolox Standards and samples, defined as the concentration at which 50% of the DPPH-radicals are scavenged.
- To determine the IC<sub>50</sub>, the concentration range of each sample should be optimized as described below.

Table 1. Amount of each solution to add to a well

	Sample	Blank 1	Blank 2
<b>Sample solution</b>	20 µL	-	-
<b>Solvent</b>	-	20 µL	20 µL
<b>Ethanol</b>	-	-	100 µL
<b>Assay Buffer</b>	80 µL	80 µL	80 µL
<b>DPPH working solution</b>	100 µL	100 µL	-

	1	2	3	4	5	6
A	Sample x 1/10 <sup>7</sup>					Blank 1
B		1/10 <sup>6</sup>				Blank 2
C		1/10 <sup>5</sup>				
D		1/10 <sup>4</sup>				
E		1/10 <sup>3</sup>				
F		1/10 <sup>2</sup>				
G		1/10				
H	Sample					

Figure 3. Example of a plate format for a preliminary experiment

\* Blank 1: coloring without antioxidant, Blank 2: sample solvent blank.

\* If the sample is highly colored, prepare a sample blank for each concentration.

### Determination of the optimum concentration range

\* Please follow the order of this protocol because it is optimised for an antioxidant assay.

1. Prepare four or more concentrations for each sample with a 10-fold dilution from the highest concentration.
2. Add 20 µL of each sample prepared in step 1 to the appropriate wells.
3. Add 20 µL of solvent used for sample extraction or dilution to the wells of Blank 1 and Blank 2.
- \* To avoid any concentration change because of volatilization, move to step 4 immediately.

4. Add 80 µL of Assay Buffer to each well.

5. Add 100 µL of ethanol to the wells of Blank 2 and mix well by pipetting.

6. Add 100 µL of DPPH working solution to the wells of the samples and Blank 1, and mix well by pipetting.

\* Because the reaction starts immediately after the addition of the DPPH working solution, it should be dispensed with a multichannel pipette to minimize any time lag in pipetting.

\* DPPH working solution is also used for the antioxidant capacity assays described later.

To avoid any concentration change because of volatilization, dispense only the required volume of DPPH working solution into a reservoir.

7. Incubate the microplate at 25°C for 30 min in the dark.

8. Measure the absorbance at 517 nm using a microplate reader.

\* If the 517 nm filter is not available, measure the absorbance at 500 nm or greater (as close as possible to 517 nm).

9. Calculate the inhibition ratio of the samples from the following equation:

$$\text{Inhibition ratio of sample (\%)} = (A_{CS} - A_S) / A_{CS} \times 100$$

A<sub>CS</sub>: Blank 1 - Blank 2

A<sub>S</sub>: Absorbance of samples - Blank 2 or sample blank (in the case where the sample is highly colored)

10. Plot the inhibition ratio (y) against the sample concentration (x) and draw a regression line (y=ax+b).

11. Determine the optimum concentration range encompassing the 50% scavenging concentration for DPPH-radicals from the regression line drawn in step 10.

Determination of the optimum concentration range

The optimum concentration range for gallic acid, a type of antioxidant, was determined as follows.

- The regression line for the concentration range of 1,000 – 0.001 µg/mL of gallic acid was plotted.
- The regression line indicated that the range between 10 and 100 µg/mL of gallic acid encompassed a 50% scavenging concentration. The IC<sub>50</sub> of gallic acid was calculated from the replotted regression line prepared from 5 to 100 µg/mL of gallic acid.

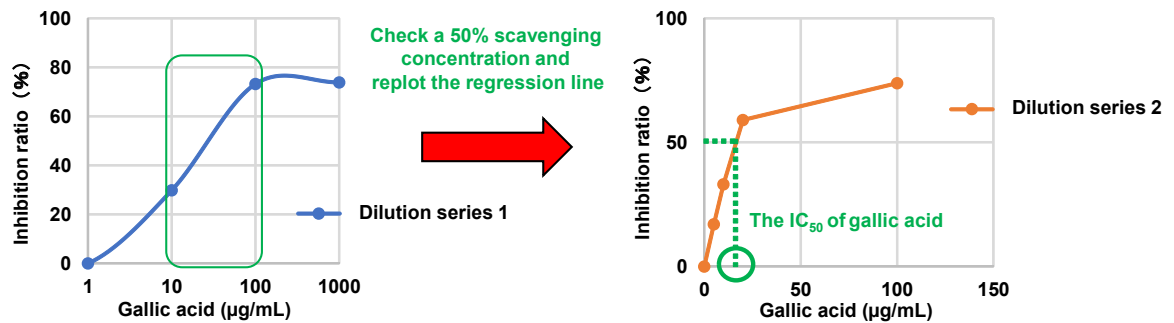


Figure 4. Example of an IC<sub>50</sub> determination for gallic acid  
 Dilution series 1: 0.001, 0.01, 0.1, 1, 10, 100, and 1000 µg/mL  
 Dilution series 2: 5, 10, 20, and 100 µg/mL

Table 2. Amount of solutions to be added

	Sample	Trolox standard	Blank 1	Blank 2	Blank 3
Sample solution	20 µL	-	-	-	-
Solvent	-	-	20 µL	20 µL	-
Ethanol	-	-	-	100 µL	120 µL
Trolox Standard solution	-	20 µL	-	-	-
Assay Buffer	80 µL	80 µL	80 µL	80 µL	80 µL
DPPH working solution	100 µL	100 µL	100 µL	-	-

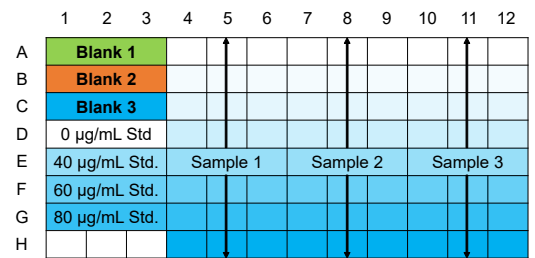


Figure 5. Example of a typical plate format

\* Blank 1: coloring without antioxidant, Blank 2: sample solvent blank, Blank 3: ethanol blank

\* If the sample is highly colored, prepare a sample blank for each concentration

(1) Measurement of the DPPH-radical scavenging ratio of Trolox and unknown samples

\* Please follow the order of this protocol because it is optimised for an antioxidant assay.

1. Add 20 µL of 0, 40, 60, and 80 µg/mL of Trolox Standard solution to each well.
  2. Add 20 µL of the sample solution at four or more concentrations (using the optimum concentration range that was determined from a preliminary experiment), to each well.
  3. Add 20 µL of ethanol to the wells of Blank 3 and add 20 µL of the solvent that was used for sample dilution to the wells of Blank 1 and Blank 2.
- \* To avoid any concentration change because of volatilization, move to step 4 immediately.
4. Add 80 µL of Assay Buffer to each well.
  5. Add 100 µL of ethanol to the wells of Blank 2 and Blank 3 and mix the wells by pipetting.
  6. Add 100 µL of DPPH working solution to the wells of Trolox, samples and Blank 1, and mix well by pipetting.
  7. Incubate the microplate at 25°C for 30 minutes in the dark.
  8. Measure the absorbance at 517 nm with a microplate reader.
- \* If a 517 nm filter is not available, measure the absorbance at 500 nm or greater (as close as possible to 517 nm).
9. Calculate the inhibition ratio of the samples from the following equation:

$$\text{Inhibition ratio of Trolox (\%)} = (A_c - A_R)/A_c \times 100$$

A<sub>c</sub>: Absorbance of 0 µg/mL Trolox Standard solution - Blank 3  
 A<sub>R</sub>: Absorbance of 40 to 80 µg/mL Trolox Standard solution - Blank 3

$$\text{Inhibition ratio of sample (\%)} = (A_{CS} - A_S)/A_{CS} \times 100$$

A<sub>CS</sub>: Blank 1 - Blank 2  
 A<sub>S</sub>: Absorbance of samples - Blank 2 or sample blank (in case the sample is highly colored)

10. Plot the inhibition ratio (y) against the sample concentration (x) and draw a regression line (y=ax+b).

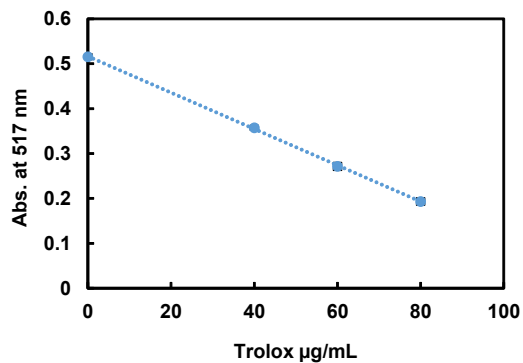


Figure 6. Absorbance changes of DPPH after Trolox treatment

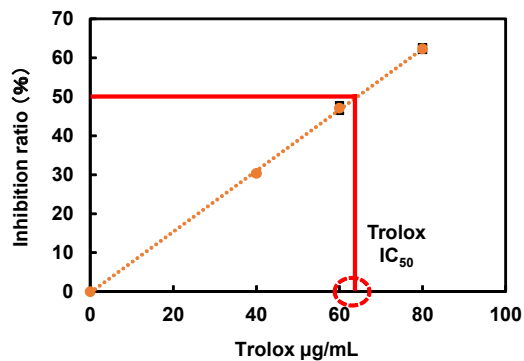


Figure 7. Inhibition ratio changes after Trolox treatment

## (2) Calculation of the Trolox equivalent antioxidant capacity (TEAC)

1. Calculate TEAC from the following equation:

$$\text{TEAC} = \text{IC}_{50}(\text{Trolox}) / \text{IC}_{50}(\text{sample})$$

1) T. Shimamura et al., *Anal. Sci.*, **2014**, 30, 717-721.

This product was commercialized under the advisory of Dr. Tomoko Shimamura (Faculty of Agriculture and Marine Science, Kochi University)

## Reference

If you need more information, please contact Dojindo technical service.

Dojindo Laboratories  
2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto  
861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525  
E-mail: info@dojindo.co.jp Web: www.dojindo.co.jp

Dojindo Molecular Technologies, Inc.  
Tel: +1-301-987-2667 Web: <http://www.dojindo.com/>  
Dojindo EU GmbH  
Tel: +49-89-3540-4805 Web: <http://www.dojindo.eu.com/>  
Dojindo China Co., Ltd  
Tel: +86-21-6427-2302 Web: <http://www.dojindo.cn/>