



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# Lipid Peroxidation (LPO) Assay

(Cat. # BAQ067)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

INTRODUCTION ..... 3

ITEM(S) SUPPLIED ..... 3

STORAGE CONDITIONS ..... 4

ADDITIONAL ITEMS REQUIRED ..... 4

SAMPLE PREPARATION ..... 4

    TISSUES AND CELLS ..... 4

    PLASMA SAMPLES..... 4

PREPARATION BEFORE USE ..... 4

    LPO REAGENT A SOLUTION..... 4

    STANDARD PREPARATION ..... 4

PROTOCOL ..... 5

DATA ANALYSIS..... 5

RELATED PRODUCTS ..... 6

## INTRODUCTION

Lipid peroxidation is a well-known example of oxidative damage in cell membranes, lipoproteins, and other lipid-containing structures. Peroxidative modification of unsaturated phospholipids, glycolipids, and cholesterol can occur in reactions triggered by i) free radical species such as oxyl radicals, peroxy radicals, and hydroxyl radicals derived from iron-mediated reduction of hydrogen peroxide or ii) non-radical species such as singlet oxygen, ozone, and peroxy nitrite generated by the reaction of superoxide with nitric oxide.

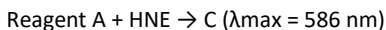
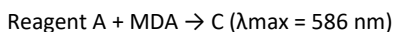
Malondialdehyde (MDA) and 4-hydroxyalkenals are important toxic byproducts of lipid peroxidation. The measurement of the amounts of such aldehydes has been widely used as an index of lipid peroxidation in vitro and in vivo.

4-Hydroxynonenal (4-HNE) is produced as a major product of the peroxidative decomposition of  $\omega$ -6 polyunsaturated fatty acids (PUFA) and possesses cytotoxic, hepatotoxic, mutagenic, and genotoxic properties. Increased levels of HNE were found in plasma and various organs under oxidative stress conditions.

MDA is in many instances the most abundant individual aldehyde resulting from lipid peroxidation. In vitro MDA can alter proteins, DNA, RNA, and many other biomolecules.

G-Biosciences' LPO assay kit measures MDA and HNE concentrations as an index of lipid peroxidation. Reactions between indoles and aldehydes (MDA and HNE) are initiated by acid-catalyzed attack at the 3-position of the indole ring to give a diindolylalkane (chromophore) with maximum absorbance in the region of 580-620 nm.

In our assay an indol (Reagent A) reacts quickly with MDA and HNE in acidic medium, yielding a chromophore (C) with a high molar extinction coefficient at its maximal absorption wavelength of 586 nm.



## ITEM(S) SUPPLIED

Description	200 tests (96 well plate)
LPO Solvent	1 vial
LPO Reagent A	3 vials (powder)
LPO Reagent B	5 vials
LPO Standard	1 vial

## STORAGE CONDITIONS

This kit is shipped at ambient temperature. Store the LPO standard at 4°C upon arrival. All other reagents can be stored at room temperature. If stored and used as directed this kit is stable for 12 months.

## ADDITIONAL ITEMS REQUIRED

- Spectrophotometer microplate reader that can measure at 580-620 nm
- 96 well microtiter plate for microplate assay.
- 1.5ml Tubes
- 20mM Tris buffer, pH 7.4

## SAMPLE PREPARATION

### *Tissues and cells*

1. Homogenize at 4°C in 20 mM Tris buffer (pH 7.4).
2. Centrifuge for 10 min at 2,000 x g and 4°C.

### *Plasma samples*

Plasma sample does not need prior sample preparation. Use directly in the assay.

## PREPARATION BEFORE USE

### *LPO Reagent A Solution*

Add 25 mL of LPO Solvent to each vial of LPO Reagent A and mix well. Once prepared, the mixture must be used immediately.

### *Standard preparation*

Prepare calibration curve in 1.5 mL tubes as shown below.

Sample	Standard [ $\mu\text{L}$ ]	H <sub>2</sub> O Diluent [ $\mu\text{L}$ ]	Standard [ $\mu\text{M}$ ]
S1 (Blank)	0	1000	0
S2	5	995	5
S3	10	990	10
S4	20	980	20
S5	30	970	30
S6	40	960	40
S7	50	950	50
S8	60	940	60

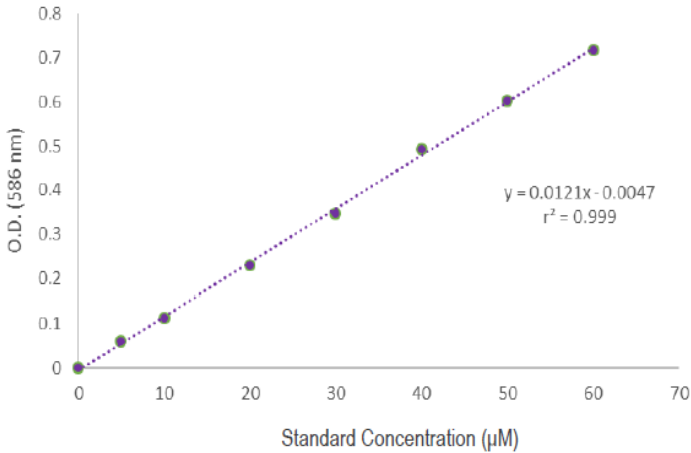
## PROTOCOL

1. Add 100  $\mu\text{L}$  of samples or standard to 325  $\mu\text{L}$  of previously prepared LPO Reagent A solution
2. Add 75  $\mu\text{L}$  of LPO Reagent B to the mixture and mix thoroughly.
3. Incubate the mixture for 40 minutes at 40°C.
4. If the mixture is cloudy, centrifuge it at 5,000 x g, for five minutes at room temperature.
5. Transfer 200  $\mu\text{L}$  of the supernatant from each tube into a 96-well plate.
6. Measure the absorbance at 586 nm using a 96-well plate reader.

## DATA ANALYSIS

Subtract the average absorbance of the Standard 1 (Blank) from itself and all other standards and samples to obtain the corrected absorbance.

Plot the corrected absorbance of standards as a function of their final concentrations.



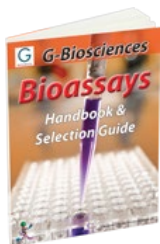
Typical standard curve for LPO assay

Calculate the aldehydes values (MDA + HNE) of the samples using the equation obtained from the linear regression of the standard curve replacing the corrected absorbance values for each sample.

$$\text{MDA + HNE } (\mu\text{M}) = (\text{sample absorbance} - y) / \text{slope}$$

## RELATED PRODUCTS

Download our Bioassays Handbook.



<http://info2.gbiosciences.com/complete-bioassay-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.



[www.GBiosciences.com](http://www.GBiosciences.com)