



Free Fatty Acid Assay Kit

Catalog Number KA1667

100 assays

Version: 05

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Intended Use	3
Background	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	4
Precautions for Use	4
Assay Protocol	5
Reagent Preparation	5
Assay Procedure	5
Data Analysis.....	7
Calculation of Results.....	7
Resources	8
References	8

Introduction

Intended Use

Applications:

- ✓ Assays: free fatty acids in biological samples such as serum, plasma, urine, saliva, milk, cell cultures and in food, agriculture products.
- ✓ Drug Discovery/Pharmacology: effects of drugs on free fatty acid metabolism.

Features

- ✓ Sensitive. Use 10 μ L samples. Linear detection range: colorimetric assay 7 - 1000 μ M, fluorimetric assay 7 - 100 μ M fatty acid.
- ✓ Convenient. Room temperature “mix-and-read” procedure can be readily automated for high-throughput assay of thousands of samples per day.

Background

Fatty acids are aliphatic monocarboxylic acids that are ubiquitously found in animal or vegetable fat, oil and wax. Fatty Acids play important roles in cellular synthesis, energy metabolism and are implicated in diverse disorders such as diabetes mellitus, sudden infant death syndrome and Reye Syndrome. Free Fatty Acid Assay Kit provides a simple, one-step and high-throughput assay for measuring free fatty acids. In this assay, free fatty acids are enzymatically converted to acyl-CoA and subsequently to H_2O_2 . The resulting H_2O_2 reacts with a specific dye to form a pink colored product. The optical density at 570nm or fluorescence intensity (530/585 nm) is directly proportional to the free fatty acid concentration in the sample.

General Information

Materials Supplied

List of component

Component	Amount
Assay Buffer	20 mL
Dye Reagent	120 µL
Enzyme A (Dried)	1 vial
Enzyme B	120 µL
CoSubstrate	120 µL
Standard	1 mL 1 mM palmitic acid

Storage Instruction

Store all components at -20°C. Shelf life of six months after receipt.

Materials Required but Not Supplied

- ✓ Pipetting devices
- ✓ Centrifuge tubes
- ✓ Clear flat-bottom uncoated 96-well plates
- ✓ Optical density plate reader
- ✓ black flat-bottom uncoated 96-well plates
- ✓ Fluorescence plate reader
- ✓ For milk and solid samples, 0.45 µm PTFE syringe filter and 5% isopropanol, 5% Triton X-100 solution.

Precautions for Use

Reagents are for research use only. Normal precautions or laboratory reagents should be exercised while using the reagents.

Assay Protocol

Reagent Preparation

Reconstitute Enzyme A by adding 120µl dH₂O to the Enzyme A tube. Make sure Enzyme A is fully dissolved by pipetting up and down and incubate at RT for 15 min. Store reconstituted Enzyme A at -20°C and use within 2 months.

Assay Procedure

✓ Colorimetric assay

Liquid samples such as serum and plasma can be assayed directly. Milk and solid samples can be homogenized in 5% isopropanol and 5% Triton X-100 in water, followed by filtration through a 0.45µm PTFE syringe filter.

Note: SH-containing reagents (e.g. β-mercaptoethanol, dithiothreitol, > 5 µM), sodium azide, EDTA, and sodium dodecyl sulfate are known to interfere in this assay and should be avoided in sample preparation.

1. Equilibrate all components to room temperature. Briefly centrifuge the tubes before opening. Keep thawed tubes on ice during assay. Important: the thawed Standard solution should be clear and colorless. If the Standard is turbid, bring it to 37°C and gently swirl the tube (do not vortex) until the solution is clear.
2. Standards: Dilute standard in Assay Buffer as follows.

No	1000 µM STD + Buffer	Vol (µL)	Palmitic Acid (µM)
1	100µL + 0µL	100	1000
2	60µL + 40µL	100	600
3	30µL + 70µL	100	300
4	0µL + 100µL	100	0

Transfer 10 µL diluted standards into separate wells of a clear flat-bottom 96-well plate.

Samples: transfer 10 µL of each sample into separate wells of the plate.

3. Color reaction. Prepare enough Working Reagent by mixing, for each well, 90 µL Assay Buffer, 1 µL Enzyme A, 1 µL Enzyme B, 1 µL CoSubstrate and 1 µL Dye Reagent. Add 90 µL Working Reagent to each well. Tap plate to mix. Incubate 30 min at room temperature.
4. Read optical density at 570 nm (550-585 nm).

✓ Fluorimetric assay

The fluorimetric assay procedure is similar to the colorimetric procedure except that (1) 0, 30, 60 and 100 µM Standards and (2) a black 96-well plate are used. Read fluorescence intensity at $\lambda_{\text{ex}} = 530 \text{ nm}$ and $\lambda_{\text{em}} = 585 \text{ nm}$.

Note: if the calculated free fatty acid concentration of a sample is higher than 1000 µM in the Colorimetric Assay or 100 µM in the Fluorimetric Assay, dilute sample in Assay Buffer and repeat the assay. Multiply

result by the dilution factor n .

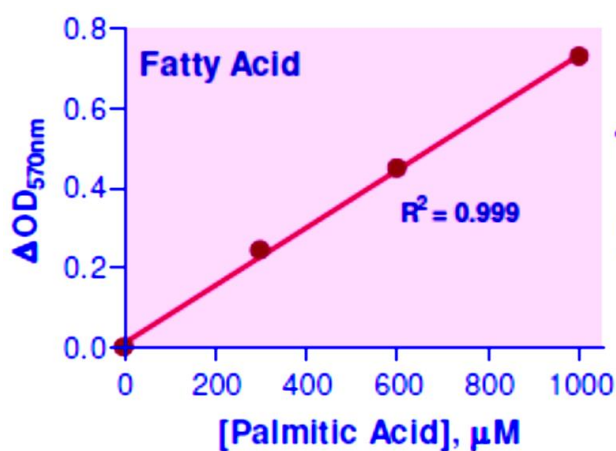
Data Analysis

Calculation of Results

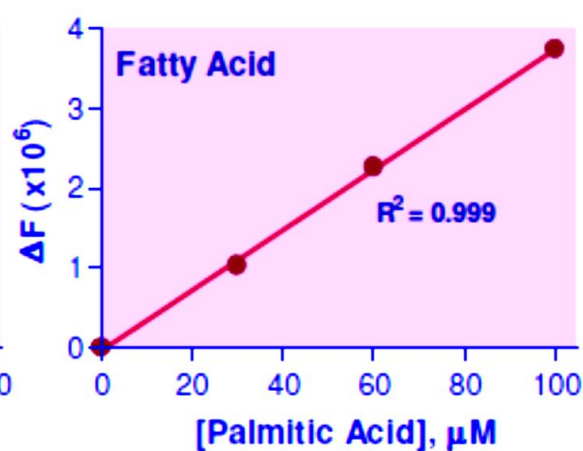
Subtract blank value (#4) from the standard values and plot the ΔOD or ΔF against standard concentrations. Determine the slope and calculate the fatty acid concentration of Sample,

$$[\text{Free Fatty Acid}] = \frac{R_{\text{SAMPLE}} - R_{\text{BLANK}}}{\text{Slope}(\mu\text{M}^{-1})} \times n(\mu\text{M})$$

R_{SAMPLE} and R_{BLANK} are optical density or fluorescence intensity readings of the Sample and Buffer Blank, respectively. n is the sample dilution factor.



96-well colorimetric assay



96-well fluorimetric assay

Resources

References

- ✓ Hwang I et al (2012) Catalase deficiency accelerates diabetic renal injury through peroxisomal dysfunction. *Diabetes* 61(3):728-738.
- ✓ Jelinek D et al (2012) The Niemann-Pick C1 gene is downregulated in livers of C57BL/6J mice by dietary fatty acids, but not dietary cholesterol, through feedback inhibition of the SREBP pathway. *J Nutr* 142(11):1935-1942.
- ✓ Seo, CW et al (2011). Antihyperlipidemic and body fat-lowering effects of silk proteins with different fibroin/sericin compositions in mice fed with high fat diet. *J Agric Food Chem* 59(8):4192-4197.