
Product Manual

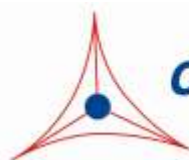
Free Fatty Acid Assay Kit (Colorimetric)

Catalog Number

STA-618

100 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.

Creating Solutions for Life Science Research

Introduction

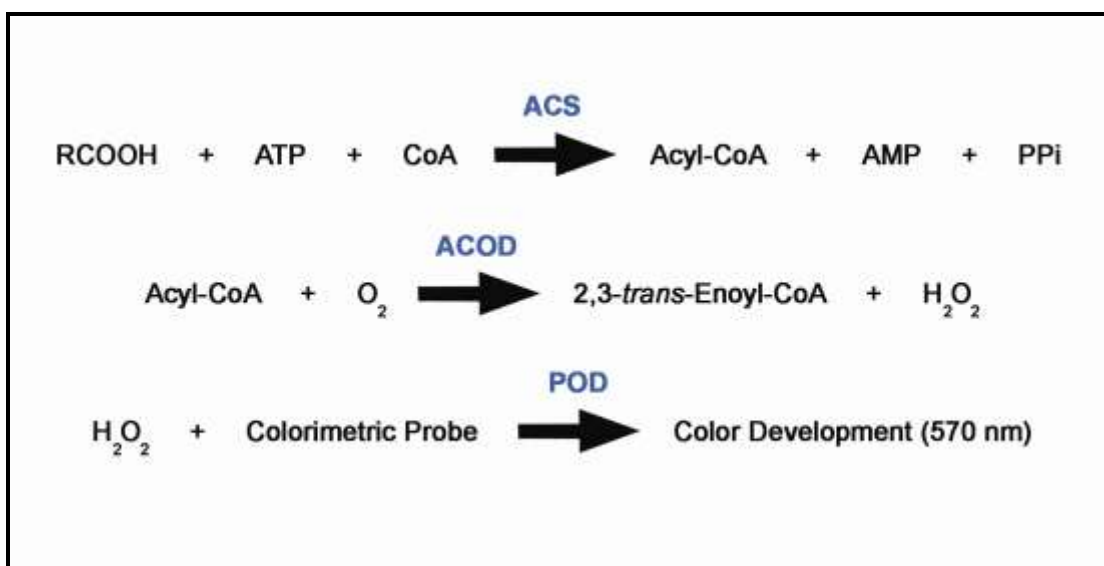
Triglycerides (TAG) are a type of lipid in the blood, serving as an energy source and playing a key role in metabolism. Triglycerides are the digestive end product of breaking down dietary fats. Any extra carbohydrates and fats that are not immediately used are chemically converted into triglycerides. In the intestines, secreted enzyme lipases hydrolyze the triglyceride ester bond, yielding glycerol and free fatty acids (FFA) in a process called lipolysis. Additionally, hormones induce and regulate lipase activity in adipose tissue, resulting in changes to blood FFA levels. Free fatty acids then bind plasma albumin for circulation in the body, serving as a readily absorbed energy source for muscle, brain and other organ tissues. Measurement of free fatty acids has become useful in monitoring and diagnosis of several diseases and metabolic disorders (e.g. obesity, insulin resistance, diabetes, cancer).

Cell Biolabs' Free Fatty Acid Assay Kit measures non-esterified fatty acids (NEFA) in serum and plasma by a coupled enzymatic reaction system (ACS-ACOD Method). First, Acyl CoA Synthetase (ACS) catalyzes fatty acid acylation of coenzyme A. Next, the acyl-CoA product is oxidized by Acyl CoA Oxidase (ACOD), producing hydrogen peroxide which reacts with the kit's Colorimetric Probe (absorbance maxima of 570 nm).

The Free Fatty Acid Assay Kit is a simple, colorimetric assay that quantitatively measures the free fatty acid concentration (non-esterified) in various samples using a 96-well microtiter plate format. Each kit provides sufficient reagents to perform up to 100 assays, including blanks, standards and unknown samples. The kit contains a palmitic acid standard and has a detection sensitivity limit of ~15 μM .

Note: This kit is not suitable for urine or heparin-containing samples. Fatty acids (C8 and longer) can be quantified with this kit.

Assay Principle



Related Products

1. STA-241: Human Low Density Lipoprotein
2. STA-243: Human High Density Lipoprotein
3. STA-369: OxiSelect™ Human Oxidized LDL ELISA Kit (MDA-LDL Quantitation)
4. STA-375: Uric Acid/Uricase Assay Kit
5. STA-378: Creatinine Assay Kit
6. STA-390: Total Cholesterol Assay Kit
7. STA-391: HDL and LDL/VLDL Cholesterol Assay Kit
8. STA-394: HDL Cholesterol Assay Kit
9. STA-396: Serum Triglyceride Quantification Kit (Colorimetric)
10. STA-397: Serum Triglyceride Quantification Kit (Fluorometric)
11. STA-398: Free Glycerol Assay Kit (Colorimetric)
12. STA-399: Free Glycerol Assay Kit (Fluorometric)

Kit Components

1. FFA Standard (Part No. 261801): One 100 µL vial of 50 mM palmitic acid in ethanol.
2. 20X Assay Buffer (Part No. 261802): One 1.5 mL vial.
3. 5X Enzyme Mixture A (Part No. 261803): Four 1 mL vials (containing ACS, Ascorbate Oxidase, and necessary cofactors).
4. 5X Enzyme Mixture B (Part No. 261804): Four 0.5 mL vials (containing ACOD).
5. NEM Reagent (Part No. 261805): One 150 µL amber vial.
6. Colorimetric Probe (Part No. 261806): One 110 µL amber vial.

Materials Not Supplied

1. 96-well microtiter plate
2. 10 µL to 1000 µL adjustable single channel micropipettes with disposable tips
3. 50 µL to 300 µL adjustable multichannel micropipette with disposable tips
4. Multichannel micropipette reservoir
5. Microplate reader capable of reading at 570 nm

Storage

Store the entire kit at -80°C. Avoid multiple freeze/thaws by aliquoting. The NEM Reagent and Colorimetric Probe are light sensitive and should be maintained in amber tubes.

Preparation of Reagents

- FFA Standard: Thaw at 37°C for 10 minutes. Once homogeneous and mixed well, maintain the standard at room temperature during assay preparation. The solution is stable for 1 week at room temperature. For longer term storage, the solution should be aliquoted and frozen at -80°C to avoid multiple freeze/thaws.
- 1X Assay Buffer: 20X Assay Buffer should be thawed/maintained at 4°C during assay preparation. Dilute the 20X Assay Buffer with deionized water. Stir to homogeneity. The 1X solution is stable for 1 week at 4°C. For longer term storage, any unused 20X stock material should be aliquoted and frozen at -80°C to avoid multiple freeze/thaws.
- 1X Enzyme Mixture (A or B): Each 5X Enzyme Mixture should be thawed/maintained at 4°C during assay preparation. Dilute the 5X Enzyme Mixture with cold, deionized water. Stir to homogeneity. The 1X solution is stable for 1 week at 4°C. For longer term storage, any unused 5X stock material should be aliquoted and frozen at -80°C to avoid multiple freeze/thaws.

Note: These components are provided in multiple tubes to minimize multiple freeze/thaws.

- NEM Reagent: Thaw and maintain at 4°C during assay preparation. The solution is stable for 1 week at 4°C. For longer term storage, the solution should be aliquoted and frozen at -80°C to avoid multiple freeze/thaws.
- Colorimetric Probe: Thaw and maintain at room temperature during assay preparation. Any unused material should be aliquoted and frozen at -80°C to avoid multiple freeze/thaws.

Preparation of FFA Standard

Heat the FFA Standard at 37°C for 10 minutes. Mix well by vortexing to ensure the solution is homogeneous and free of palmitic acid crystals. Freshly prepare a dilution series of standard in the concentration range of 0 μM – 500 μM by diluting the standard stock solution (provided at 50 mM) in 1X Assay Buffer (see Table 1). FFA diluted solutions and standards should be prepared fresh, vortexed well and used immediately.

Standard Tubes	50 mM FFA Standard (μL)	1X Assay Buffer (μL)	Final FFA Standard (μM)
1	10	990	500
2	500 of Tube #1	500	250
3	500 of Tube #2	500	125
4	500 of Tube #3	500	62.5
5	500 of Tube #4	500	31.25
6	500 of Tube #5	500	15.63
7	500 of Tube #6	500	7.81
8	0	500	0

Table 1. Preparation of Free Fatty Acid Standards

Preparation of Samples

- Urine: This kit is not recommended for urine samples.
- Plasma: Collect blood with an anticoagulant such as citrate, EDTA or oxalate and mix by inversion. Centrifuge the blood at 1000 x g at 4°C for 10 minutes. Collect plasma supernatant without disturbing the white buffy layer. Sample should be tested immediately or frozen at -80°C for storage. Plasma must be diluted before assaying (1:2 to 1:20 in 1X Assay Buffer). Normal FFA levels in human plasma are typically 150-450 µM.

Note: Heparin is known to interfere with the assay. Heparin-containing samples including heparinized plasma should be avoided.

- Serum: Collect blood in a tube with no anticoagulant. Allow the blood to clot at room temperature for 30 minutes. Centrifuge at 2500 x g for 20 minutes. Remove the yellow serum supernatant without disturbing the white buffy layer. Samples should be tested immediately or frozen at -80°C for storage. Serum must be diluted before assaying (1:2 to 1:20 in 1X Assay Buffer). Normal FFA levels in human serum are typically 100-700 µM.

Assay Protocol

Each FFA standard and sample should be assayed in duplicate or triplicate. A freshly prepared standard curve should be used each time the assay is performed.

1. Add 10 µL of each FFA standard, sample and blank to the 96-well microtiter plate.
2. Add 200 µL of 1X Enzyme Mixture A (see Preparation of Reagents) to each well.
3. Cover the plate wells to protect the reaction from light.
4. Incubate at 37°C for 30 minutes.
5. During the step 4 incubation, separately prepare the desired volume of Detection Enzyme Mixture according to Table 2 below, based on the number of tests to be performed. Maintaining all components and mixtures at 4°C throughout this step, add components in the following sequence:
 - a. In a tube, add the appropriate volume of 1X Enzyme Mixture B (see Preparation of Reagents).
 - b. To the 1X Enzyme Mixture B add the corresponding volume of NEM Reagent. Mix well.
 - c. Finally, add the corresponding volume of Colorimetric Probe. Mix well and use immediately.

Note: Detection Enzyme Mixture will appear slightly pink in color. This is normal background and should be subtracted from all absorbance values.

1X Enzyme Mixture B (mL)	NEM Reagent (µL)	Colorimetric Probe (µL)	Total Volume of Detection Enzyme Mixture (mL)	# of Tests in 96-well Plate (100 µL/test)
10	100	100	10.2	100
5	50	50	5.1	50
2.5	25	25	2.55	25

Table 2. Preparation of Detection Enzyme Mixture

- Transfer 100 μL of the above Detection Enzyme Mixture to each well (from step 4).
- Cover the plate wells to protect the reaction from light.
- Incubate at 37°C for 10 minutes.
- Read absorbance in the 540-570 nm range on a microplate reader.
- Calculate the concentration of free fatty acid within samples by comparing the sample absorbance to the standard curve. Negative controls (without FFA) should be subtracted.

Example of Results

The following figures demonstrate typical Free Fatty Acid Assay Kit results. One should use the data below for reference only. This data should not be used to interpret actual results.

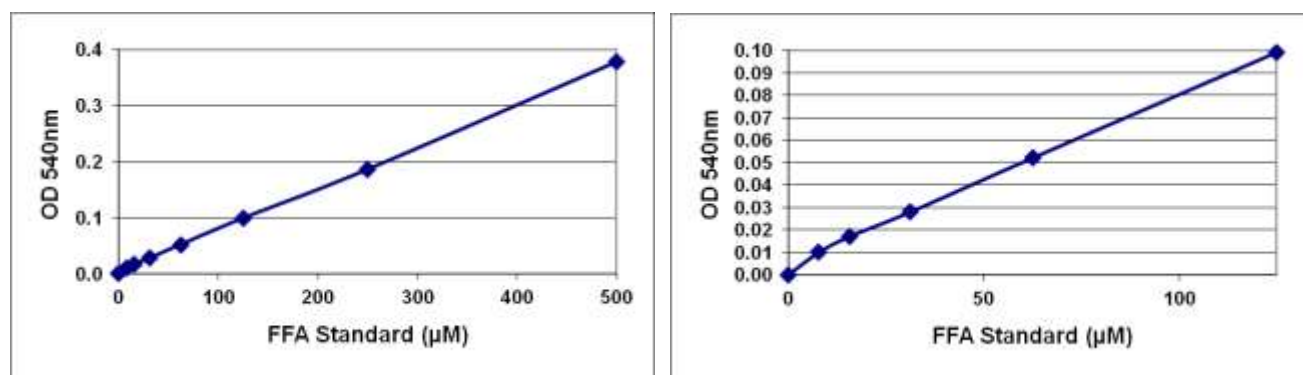


Figure 1: Free Fatty Acid Assay Standard Curve. FFA standard curve was performed according to the Assay Protocol. Background has been subtracted.

References

- Boden, G., Cheung, P., Stein, T.P. et al. (2002) *Am J. Physiol.* **283**, E12-E-19.
- Steinberg, H.O., Tarshoby, M., Monestel, R. et al. (1997) *J. Clin. Invest.* **100**, 1230-1239.
- Boden, G., Lebed, B. Schatz, M. et al. (2001) *Diabetes* **50**, 1612-1617.
- Boden, G., She, P., Mozzoli, M. et al. (2005) *Diabetes* **54**, 917-927.
- Kelley, D.E., Mokan, M., Simoneau, J.A., and Mandarino, L.J. (1993) *J. Clin. Invest.* **92**, 91-98.
- Lewis, G.F., Uffelman, K.D., Szeto, L.W., Weller, B., and Steiner, G. (1995) *J. Clin. Invest.* **95**, 158-166.

Recent Product Citations

- Devarshi, P.P. et al. (2021). A single bout of cycling exercise induces nucleosome repositioning in the skeletal muscle of lean and overweight/obese individuals. *Diabetes Obes Metab.* doi: 10.1111/dom.14541.
- Valladolid-Acebes, I. et al. (2021). Lowering apolipoprotein CIII protects against high-fat diet-induced metabolic derangements. *Sci Adv.* **7**(11):eabc2931. doi: 10.1126/sciadv.abc2931.
- Casey, C.A. et al. (2021). Lipid droplet membrane proteome remodeling parallels ethanol-induced hepatic steatosis and its resolution. *J Lipid Res.* doi: 10.1016/j.jlr.2021.100049.
- Zhang, J. et al. (2020). ADORA1-driven brain-sympathetic neuro-adipose connections control body weight and adipose lipid metabolism. *Mol Psychiatry.* doi: 10.1038/s41380-020-00908-y.

5. Kale, M. et al. (2020). Modulates Anxiety and Depression-like Behaviour in Diabetic Insulin-Resistant Rats. *Brain Res.* doi: 10.1016/j.brainres.2020.147045.
6. Miller, C.N. et al. (2019). Fetal growth outcomes following peri-implantation exposure of Long-Evans rats to noise and ozone differ by sex. *Biol Sex Differ.* **10**(1):54. doi: 10.1186/s13293-019-0270-6.
7. Henriquez, A.R. et al. (2019). Exacerbation of ozone-induced pulmonary and systemic effects by β 2-adrenergic and/or glucocorticoid receptor agonist/s. *Sci Rep.* **9**(1):17925. doi: 10.1038/s41598-019-54269-w.
8. Martinez, N. et al. (2019). mTORC2/Akt activation in adipocytes is required for adipose tissue inflammation in tuberculosis. *EBioMedicine.* pii: S2352-3964(19)30433-5. doi: 10.1016/j.ebiom.2019.06.052.
9. Matsue, M. et al. (2019). Measuring the Antimicrobial Activity of Lauric Acid against Various Bacteria in Human Gut Microbiota Using a New Method. *Cell Transplant.* doi: 10.1177/0963689719881366.
10. Tillman, M.C. et al. (2019). Structural characterization of life-extending *Caenorhabditis elegans* Lipid Binding Protein 8. *Sci Rep.* **9**(1):9966. doi: 10.1038/s41598-019-46230-8.
11. Miller, C.N. et al. (2019). Ozone Exposure During Implantation Increases Serum Bioactivity in HTR-8/SVneo Trophoblasts. *Toxicol Sci.* **168**(2):535-550. doi: 10.1093/toxsci/kfz003.
12. Miller, C.N. et al. (2019). Aspirin pre-treatment modulates ozone-induced fetal growth restriction and alterations in uterine blood flow in rats. *Reprod Toxicol.* **83**:63-72. doi: 10.1016/j.reprotox.2018.12.002.
13. Allard, C. et al. (2019). Loss of Nuclear and Membrane Estrogen Receptor- α Differentially Impairs Insulin Secretion and Action in Male and Female Mice. *Diabetes.* **68**(3):490-501. doi: 10.2337/db18-0293.
14. Reijnders, D. et al. (2019). Dyslipidemia and the role of the adipose tissue in early pregnancy in the BPH/5 mouse model for preeclampsia. *Am J Physiol Regul Integr Comp Physiol.* doi: 10.1152/ajpregu.00334.2018.
15. El-Shiekh, R.A. et al. (2019). Anti-obesity effect of argel (*Solenostemma argel*) on obese rats fed a high fat diet. *J Ethnopharmacol.* **238**:111893. doi: 10.1016/j.jep.2019.111893.
16. Attia, R.T. et al. (2019). Raspberry ketone and Garcinia Cambogia rebalanced disrupted insulin resistance and leptin signaling in rats fed high fat fructose diet. *Biomed Pharmacother.* **110**:500-509. doi: 10.1016/j.biopha.2018.11.079.
17. Paris, H.L. et al. (2019). Effect of carbohydrate ingestion on central fatigue during prolonged running exercise in moderate hypoxia. *J Appl Physiol (1985).* **126**(1):141-151. doi: 10.1152/jappphysiol.00684.2018.
18. Xiao, W.C. et al. (2018). Alleviation of palmitic acid-induced endoplasmic reticulum stress by augmenter of liver regeneration through IP3R-controlled Ca²⁺ release. *J Cell Physiol.* **233**(8):6148-6157. doi: 10.1002/jcp.26463.
19. Martin, B.L. et al. (2018). Acute peat smoke inhalation sensitizes rats to the postprandial cardiometabolic effects of a high fat oral load. *Sci Total Environ.* **643**:378-391. doi: 10.1016/j.scitotenv.2018.06.089.
20. Tomic, M. et al. (2018). Lsd1 regulates skeletal muscle regeneration and directs the fate of satellite cells. *Nat Commun.* **9**(1):366. doi: 10.1038/s41467-017-02740-5.
21. Matuszek, M.A. et al. (2018). Statins Do Not Impair Whole-body Fat Oxidation During Moderate-intensity Exercise in Dyslipidemic Adults. *Exerc Med.* **2**:12. doi: 10.26644/em.2018.012.

22. Hyatt, H.W. et al. (2017). Lactation has persistent effects on a mother's metabolism and mitochondrial function. *Sci Rep.* **7**(1):17118. doi: 10.1038/s41598-017-17418-7.
23. Miller, C.N. et al. (2017). Uterine Artery Flow and Offspring Growth in Long-Evans Rats following Maternal Exposure to Ozone during Implantation. *Environ Health Perspect.* **125**(12):127005. doi: 10.1289/EHP2019.
24. Xiao, X., et al. (2017). Permethrin alters glucose metabolism in conjunction with high fat diet by potentiating insulin resistance and decreases voluntary activities in female C57BL/6J mice. *Food Chem Toxicol.* **108** (Pt A):161-170. doi: 10.1016/j.fct.2017.07.053
25. Passlack, N. et al. (2017). Impact of hyperlipidaemia on intermediary metabolism, faecal microbial metabolites and urinary characteristics of lipoprotein lipase deficient vs. normal cats. *J. Anim. Physiol. Anim. Nutr. (Berl)* doi:10.1111/jpn.12721.
26. Miller, D. B. et al. (2016). Systemic metabolic derangement, pulmonary effects, and insulin insufficiency following subchronic ozone exposure in rats. *Toxicol Appl Pharmacol.* doi: 10.1016/j.taap.2016.06.027.
27. Healy, M. E. et al. (2016). Dietary sugar intake increases liver tumor incidence in female mice. *Sci Rep.* doi:10.1038/srep22292.
28. Nicholas, D. A. et al. (2015). Identification of anti-long chain saturated fatty acid IgG antibodies in serum of patients with type 2 diabetes. *Mediators Inflamm.* doi:10.1155/2015/196297.
29. Kahouli, I. et al. (2015). In-vitro characterization of the anti-cancer activity of the probiotic bacterium *Lactobacillus fermentum* NCIMB 5221 and potential against colorectal cancer. *J Cancer Sci Ther.* **7**:224-235.
30. Ditzel, E. J. et al. (2015). Effects of arsenite exposure during fetal development on energy metabolism and susceptibility to diet-induced fatty liver disease in male mice. *Environ Health Perspect.* doi:10.1289/ehp.1409501.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
 7758 Arjons Drive
 San Diego, CA 92126
 Worldwide: +1 858-271-6500
 USA Toll-Free: 1-888-CBL-0505
 E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2013-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.