



Triglyceride Quantification Kit

Catalog Number KA0847

100 assays

Version: 06

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Background	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Assay Protocol	5
Reagent Preparation	5
Sample Preparation.....	5
Assay Procedure	6
Data Analysis.....	7
Calculation of Results.....	7

Introduction

Background

Triglycerides (TG) are the main constituent of vegetable oil, animal fat, LDL and VLDL, and play an important role as transporters of fatty acids as well as serving as an energy source. TG are broken down into fatty acids and glycerol, after which both can serve as substrates for energy producing and metabolic pathways. High blood levels of TG are implicated in atherosclerosis, heart disease and stroke as well as in pancreatitis. The Triglyceride Quantification Kit provides a sensitive, easy assay to measure TG concentration in a variety of samples. In the assay, TG are converted to free fatty acids and glycerol. The glycerol is then oxidized to generate a product which reacts with the probe to generate color (spectrophotometry at $\lambda = 570 \text{ nm}$) and fluorometric (Ex/Em = 535/587 nm). The kit can detect 2 pmol-10 nmol (or 2-10000 μM range) of triglyceride in various samples. The kit also detects monoglycerides and diglycerides.

General Information

Materials Supplied

List of component

Component	Amount
Triglyceride Assay Buffer	25 ml
Triglyceride Probe (in DMSO, anhydrous)	200 µl
Lipase	1 vial
Triglyceride Enzyme Mix (lyophilized)	1 vial
Triglyceride Standard (1 mM)	0.3 ml

Storage Instruction

Store kit at -20°C, protect from light. Warm Triglyceride Assay Buffer to room temperature before use. Briefly centrifuge all small vials prior to opening.

Assay Protocol

Reagent Preparation

- ✓ Triglyceride standard: Frozen storage may cause the triglyceride standard to separate from the aqueous phase. To re-dissolve, keep the cap tightly closed, place in a hot water bath (~80-100°C) for 1 min or until the standard looks cloudy, vortex for 30 sec, the standard should become clear. Repeat the heat and vortex one more time. The triglyceride standard is now completely in solution and ready to be used.
- ✓ Triglyceride Probe: Ready to use as supplied. Warm to by placing in a 37 °C bath for 1-5 min to thaw the DMSO solution before use. (*Note: DMSO tends to be a solid after -20 °C storage, even when left at room temperature-so need to melt for a few minutes at 37 °C*). Store at -20 °C, protect from light. Use within two months.
- ✓ Triglyceride Enzyme Mix: Dissolve in 220 µl Triglyceride Assay Buffer. Aliquot and store at -20 °C. Use within two months.
- ✓ Lipase: Dissolve in 220 µl Triglyceride Assay Buffer. Aliquot and store at -20 °C. Use within two months.

Sample Preparation

Add 2-50 µl test samples to a 96-well plate. Adjust the volume to 50 µl/well with Triglyceride Assay Buffer. We suggest using dilutions volumes of sample to ensure readings are within the standard curve range. A background control should be performed by replacing 2 µl Lipase with 2 µl Triglyceride Assay Buffer. The background should be subtracted from all readings. Endogenous compounds in the sample may interfere with the assay, so it is suggested to spike samples with a known amount of Standard (4 nmol) to ensure accurate determinations of Triglyceride in your sample.

**Notes: Serum contains 0.1-6 mM triglyceride, which can be tested directly. For tissues (~100 mg), cells (~10 million) or other non-aqueous samples, homogenize in 1 ml solution containing 5% NP-40 in water, slowly heat the samples to 80-100 °C in a water bath for 2-5 minutes or until the NP-40 becomes cloudy, then cool down to room temperature. Repeat the heating one more time to solublize all triglyceride. Centrifuge for 2 min. (top speed using a microcentrifuge) to remove any insoluble materials. Dilute 10 folds with dH₂O before the assay.*

Assay Procedure

1. Standard Curve Preparation

For the colorimetric assay, Dilute 40 μ l of the 1 mM Triglyceride into 160 μ l Triglyceride Assay Buffer, mix to generate 0.2 mM Triglyceride standard. Add 0, 10, 20, 30, 40, 50 μ l of the 0.2 mM Triglyceride Standard into a series of wells. Adjust volume to 50 μ l/well with Triglyceride Assay Buffer to generate 0, 2, 4, 6, 8, 10 nmol/well of Triglyceride Standard.

For the fluorometric assay, dilute the Triglyceride Standard to 0.02 mM with the Triglyceride Assay Buffer (Detection sensitivity is 10-100 fold higher for a fluorometric than a colorimetric assay). Follow the procedure as the colorimetric assay.

2. Lipase: Add 2 μ l of lipase to each standard and sample well. Mix and incubate 20 min at room temperature to convert triglyceride to glycerol and fatty acid.

Note: If samples contain glycerol, do a sample background control, omit the lipase to determine glycerol background only, not triglyceride.

3. Triglyceride Reaction Mix: Mix enough reagent for the number of samples and standards to be performed: For each well, prepare a total 50 μ l Reaction Mix:

	Reaction Mix
Triglyceride Assay Buffer	46 μ l
Triglyceride Probe*	2 μ l
Triglyceride Enzyme Mix	2 μ l

Add 50 μ l of the Reaction Mix to each well containing the Triglyceride Standard, samples and background control(s). Mix well. Incubate at room temperature for 30-60 min. (60 min. gives slightly better result) protect from light.

**Note: For the fluorometric assay, use 0.4 μ l/well of the Probe to decrease the background readings, therefore increase detection sensitivity.*

4. Measurement: Measure absorbance at 570 nm for colorimetric assay or fluorescence at Ex/Em = 535/590 nm for fluorometric assay in a microtiter plate reader. The reaction is stable for at least two hr.

Data Analysis

Calculation of Results

Calculations: Subtract 0 Standard reading from all readings. If sample background control reading is significant then subtract the sample background control reading from sample reading. Plot the TG Standard Curve. For unspiked samples, apply the corrected OD to the TG Standard Curve to get B nmol of TG in the sample well.

Sample TG concentration (C) = $B/V \times D$ nmol/ μ l or mM

Where: B is the amount of TG from Standard Curve (nmol)

V is the sample volume added into the reaction well (μ l)

D is the sample dilution factor

Note: For spiked samples, correct for any sample interference by subtracting the sample reading from spiked sample reading.

Triglyceride molecular weight: 885.4 g/mol