



Free beta-CG (Human) ELISA Kit

Catalog Number KA0210

96 assays

Version: 04

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Intended Use	3
Background	3
Principle of the Assay	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	4
Precautions for Use	4
Assay Protocol	6
Reagent Preparation	6
Sample Preparation	6
Assay Procedure	6
Data Analysis	8
Calculation of Results	8
Performance Characteristics	9
Resources	13
References	13
Plate Layout	15

Introduction

Intended Use

For the quantitative determination of free beta subunit of human chorionic gonadotropin (free β -hCG) concentration in human serum.

Background

Human Chorionic Gonadotropin (hCG) is a glycoprotein hormone normally produced by placenta during pregnancy. The hormone is present in blood and urine around seven to thirteen days following implantation of the fertilized ovum. Structurally intact hCG molecules consist of two non-covalently linked polypeptide subunits, the alpha and beta chain subunits. Measurement of intact hCG and of the alpha subunit of hCG appears to give similar results in blood and urine but not the levels of beta subunit. In the normal secondtrimester maternal sera, the level of intact hCG range from 20,000 mIU/mL to 50,000 mIU/mL (1 ng = 15 mIU). In contrast, the levels of either free α - or free β -hCG are on average one half of 1% of hCG levels. hCG and the free subunits appear not to be useful as serological markers for nontrophoblastic tumors; however, the absolute increase of β -hCG level in choriocarcinoma patients clearly differentiates it from normal pregnancy.

Principle of the Assay

The Free beta-CG (Human) ELISA Kit is based on the principle of a solid phase enzyme-linked immunosorbent assay.¹ The assay system utilizes a unique monoclonal antibody directed against a distinct antigenic determinant on the free β -hCG. Mouse monoclonal anti-free- β -hCG antibody is used for solid phase immobilization (on the microtiter wells). A goat anti whole hCG antibody is in the antibody-enzyme (horseradish peroxidase) conjugate solution. The test sample is allowed to react sequentially with the two antibodies, resulting in the free β -hCG molecules being sandwiched between the solid phase and enzyme-linked antibodies. After two separate 30 minute incubations at 37 °C, the wells are washed with water to remove unbound labeled antibodies. A solution of TMB Reagent is added and incubated for 20 minutes, resulting in the development of a blue color.

The color development is stopped with the addition of Stop Solution changing the color to yellow. The concentration of β -hCG is directly proportional to the color intensity of the test sample. Absorbance is measured spectrophotometrically at 450 nm.

General Information

Materials Supplied

List of component

Component	Amount
Antibody-Coated Wells: Microtiter wells coated with monoclonal anti-free- β -hCG.	96 wells.
Reference Standard Set: Contains 0, 2.5, 5, 10, 25, and 50 ng/mL of β -hCG in bovine serum with preservatives, lyophilized.	1 vial x 6
Zero Buffer: Contains tris buffer with preservatives.	13 mL
Enzyme Conjugate Reagent: Contains goat anti-whole hCG conjugated to horseradish peroxidase with preservatives.	18 mL
TMB Reagent: Contains 3, 3', 5, 5' tetramethylbenzidine (TMB) stabilized in buffer solution.	11 mL
Stop Solution (1 N HCl): Diluted hydrochloric acid.	11 mL

Storage Instruction

- ✓ Store the unopened kit at 2-8°C upon receipt and when it is not in use, until the expiration shown on the kit label. Refer to the package label for the expiration date.
- ✓ Keep microtiter plate in a sealed bag with desiccant to minimize exposure to damp air.

Materials Required but Not Supplied

- ✓ Precision pipettes: 50 μ L, 100 μ L, 150 μ L, and 1.0 mL.
- ✓ Disposable pipette tips.
- ✓ Distilled water.
- ✓ Vortex mixer or equivalent.
- ✓ Absorbent paper or paper towel.
- ✓ Graph paper.
- ✓ Microtiter plate reader: A microtiter well reader with a bandwidth of 10 nm or less and an optical density range of 0 to 3 OD or greater at 450 nm wavelength is acceptable for absorbance measurement.

Precautions for Use

- Warning and Precautions
- ✓ CAUTION: This kit contains human material. The source material used for manufacture of this component tested negative for HBsAg, HIV 1/2 and HCV by FDA-approved methods. However, no method can completely assure absence of these agents. Therefore, all human blood products, including serum samples, should be considered potentially infectious. Handling should be as defined by an

appropriate national biohazard safety guideline or regulation, where it exists.²⁻⁴

- ✓ Avoid contact with 1 N HCl. It may cause skin irritation and burns. If contact occurs, wash with copious amounts of water and seek medical attention if irritation persists.
- ✓ Do not use reagents after expiration date and do not mix or use components from kits with different lot numbers.
- ✓ Replace caps on reagents immediately. Do not switch caps.
- ✓ Do not pipette reagents by mouth.
- ✓ For in Research use.

- Limitation of procedures
- ✓ Reliable and reproducible results will be obtained when the assay procedure is carried out with a complete understanding of the package insert instructions and with adherence to good laboratory practice.
- ✓ The wash procedure is critical. Insufficient washing will result in poor precision and falsely elevated absorbance readings.
- ✓ Serum samples demonstrating gross lipemia, gross hemolysis, or turbidity should not be used with this test.
- ✓ For professional use only. The results obtained from the use of this kit should be used only as an adjunct to other procedures and information available to the physician.

Assay Protocol

Reagent Preparation

- ✓ All reagents should be brought to room temperature (18-25°C) before use.
- ✓ Reconstitute each lyophilized standard with 1.0 mL distilled water. Allow the reconstituted material to stand for at least 20 minutes and mix gently. Reconstituted standards will be stable for up to 30 days when stored sealed at 2-8°C.

Sample Preparation

- The use of serum samples is required for this test.
- Specimens should be collected using standard venipuncture techniques. Remove serum from the coagulated or packed cells within 60 minutes after collection.
- Specimens which cannot be assayed within 24 hours of collection should be frozen at –20°C or lower, and will be stable for up to six months.
- Avoid grossly hemolytic (bright red), lipemic (milky), or turbid samples (after centrifugation).
- Specimens should not be repeatedly frozen and thawed prior to testing. DO NOT store in “frost free” freezers, which may cause occasional thawing. Specimens which have been frozen, and those which are turbid and/or contain particulate matter, must be centrifuged prior to use.

Assay Procedure

Procedural Notes:

- ✓ *Pipetting Recommendations (single and multi-channel): Pipetting of all standards, samples, and controls should be completed within 3 minutes.*
 - ✓ *All standards, samples, and controls should be run in duplicate concurrently so that all conditions of testing are the same.*
 - ✓ *It is recommended that the wells be read within 15 minutes following addition of Stop Solution.*
1. Secure the desired number of coated wells in the holder.
 2. Dispense 50 µL of standards, specimens, and controls into appropriate wells.
 3. Dispense 100 µL of Zero Buffer into each well.
 4. Thoroughly mix for 30 seconds. It is very important to mix them completely.
 5. Incubate at 37°C for 30 minutes.
 6. Remove the incubation mixture by flicking plate contents into a sink.
 7. Rinse and flick the microtiter wells 5 times with distilled or deionized water. (Please do not use tap water.)
 8. Strike the wells sharply onto absorbent paper or paper towels to remove all residual water droplets.
 9. Dispense 150 µL of Enzyme Conjugate Reagent into each well. Gently mix for 10 seconds.

10. Incubate at 37°C for 30 minutes.
11. Remove the incubation mixture by flicking plate contents into a waste container.
12. Rinse and flick the microtiter wells 5 times with distilled or deionized water. (Please do not use tap water.)
13. Strike the wells sharply onto absorbent paper or paper towels to remove all residual water droplets.
14. Dispense 100 μ L of TMB Reagent into each well. Gently mix for 10 seconds.
15. Incubate at room temperature for 20 minutes.
16. Stop the reaction by adding 100 μ L of Stop Solution to each well.
17. Gently mix for 30 seconds. It is important to make sure that all the blue color changes to yellow color completely.
18. Read the optical density at 450 nm with a microtiter well reader within 15 minutes.

Data Analysis

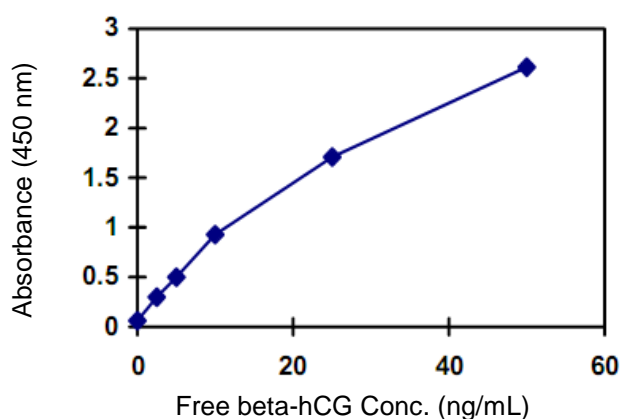
Calculation of Results

- ✓ Calculate the mean absorbance values (A_{450}) for each set of reference standards, controls, and samples.
- ✓ Construct a standard curve by plotting the mean absorbance obtained for each reference standard against its concentration in ng/mL on graph paper, with absorbance values on the vertical or Y axis, and concentrations on the horizontal or X axis.
- ✓ Use the mean absorbance values for each specimen to determine the corresponding concentration of free β -hCG in ng/mL from the standard curve.

- Example of standard curve

Results of a typical standard run with optical density readings at 450 nm shown in the Y axis against free β -hCG concentrations shown in the X axis. This standard curve is for the purpose of illustration only, and should not be used to calculate unknowns. Each user should obtain his or her own data and standard curve.

β -hCG(ng/mL)	Absorbance (450 nm)
0	0.061
2.5	0.296
5.0	0.498
10.0	0.929
25.0	1.711
50.0	2.613



Performance Characteristics

- **Quality Control**

Good laboratory practice requires that quality control specimens (controls) be run with each calibration curve to verify assay performance. To ensure proper performance, control material should be assayed repeatedly to establish mean values and acceptable ranges.

- **Expected Values**

The following information is cited from references #9, 11, 13, 14, and 15:

1. **hCG and Free β -hCG Subunit Levels in Normal Pregnancy**

A logarithmic increase in the serum concentration of hCG was observed from 5-8 weeks of gestation (2,600 ng/mL to 33,000 ng/mL) as defined by last menstrual period; thereafter, hCG values decreased. Similarly, free β -hCG levels increased rapidly to reach maximum levels (~60 ng/mL) at 8-9 weeks of pregnancy, followed by a gradual decline during the next 11-12 weeks of gestation.

At 5 weeks of gestation, the ratio of free β -hCG to intact hCG is approximately 1.0 % (w/w). Thereafter, this ratio remains remarkably constant over 22 weeks of gestation (~0.5 % w/w).

2. **hCG and Free β -hCG Subunit Levels in Gestational Choriocarcinoma**

Free α and free β -subunits and hCG levels were measured in five individual with untreated gestational choriocarcinoma. The concentrations in serum are shown in the following table:

Sample Number	hCG (ng/mL)	α -hCG (ng/mL)	β -hCG (ng/mL)
1	210,000	112	8,000
2	22,195	20	1,300
3	6,840	1	232
4	36,000	44	3,900
5	4,200	2	350

The levels of free α -hCG were low, ranging from 1-112 ng/mL, whereas hCG levels ranged from 4,200 to 210,000 ng/mL (1 ng \approx 15 mIU). In contrast, free β -hCG concentrations were found to be markedly elevated in choriocarcinoma.

- **Sensitivity**

The minimum detectable concentration of free β -hCG by this assay is estimated to be 0.25 ng/mL.

- **Precision**

- a. **Intra-Assay Precision**

Within-run precision was determined by replicate determinations of three different control sera in one assay.

Within-assay variability is shown below:

Serum Sample	1	2	3
Number of Replicates	24	24	24
Mean free β -hCG (ng/mL)	2.8	14	36
Standard Deviation	0.11	0.4	1.5
Coefficient of Variation(%)	3.7	2.6	4.3

b. Inter-Assay Precision

Between-run precision was determined by replicate measurements of three different control sera in several different assays. Between-assay variability is shown below:

Serum Sample	1	2	3
Number of Replicates	20	20	20
Mean free β -hCG (ng/mL)	3.0	17	37
Standard Deviation	0.10	0.7	0.7
Coefficient of Variation(%)	3.4	4.0	2.0

- Recovery and Linearity Studies

a. Recovery

Various serum samples of known free β -hCG levels were mixed and assayed in duplicate. The average recovery was 96.7%.

	Expected Concentration (ng/mL)	Observed Concentration (ng/mL)	% Recovery
1	41.22	39.68	96.3
2	40.61	37.73	92.9
3	15.07	15.58	103.4
4	17.44	17.06	97.8
5	3.389	3.258	96.1
6	3.352	2.986	89.1
7	0.493	0.478	97.1
8	0.436	0.441	101.1
	Average Recovery= 96.7%		

b. Linearity

Four samples were serially diluted with the zero standard in a linearity study. The average recovery was 111.3%.

#	Dilution	Expected Conc. (ng/mL)	Observed Conc. (ng/mL)	% Recovery
1.	Undiluted	-----	34.54	-----
	1:2	17.27	19.18	111.1
	1:4	8.634	8.822	102.2
	1:8	4.317	4.701	108.9
	1:16	2.159	2.398	111.1
	Mean = 108.3%			
2.	Undiluted	-----	47.81	-----
	1:2	23.91	23.98	100.3
	1:4	11.95	13.69	114.5
	1:8	5.977	7.538	126.1
	1:16	2.988	3.872	129.6
	1:32	1.494	1.921	128.6
	Mean = 117.6%			
3.	Undiluted	-----	42.09	-----
	1:2	21.05	22.70	107.9
	1:4	10.52	10.50	99.8
	1:8	5.262	5.519	104.9
	1:16	2.631	2.930	111.4
	1:32	1.315	1.622	123.3
	Mean = 106.0%			
4.	Undiluted	-----	35.48	-----
	1:2	17.74	20.27	114.3
	1:4	8.869	9.424	106.3
	1:8	4.435	5.200	117.3
	1:16	2.217	2.554	115.2
	1:32	1.109	1.367	123.3
	Mean = 113.3%			

- Specificity

The following substances were tested for cross-reactivity:

Analytes and Concentration	% Cross Reactivity
Intact hCG (100,000 mIU/mL)	0.5 %
Beta-hCG	100.0 %
Alpha-hCG (500 ng/mL)	0.0 %
LH (500 mIU/mL)	0.0 %
TSH (500 μ IU/mL)	0.0 %
FSH (500 mIU/mL)	0.0 %

- Standardization

For intact hCG, 1 ng is approximately equivalent to 15 mIU (WHO, 1st IRP 75/537). For free β -hCG subunit, since there is no WHO standardization, we tested the free β -hCG against Abnova's hCG ELISA kit, and found 1 ng of free β -hCG equals to 0.1 mIU in terms of hCG immunological activity.

Resources

References

1. Engall, E., Enzyme immunoassay ELISA and EMIT. In: Van Vunakis, H. and Langone, J.J. (eds.), *Methods in Enzymology*, Academic Press, New York, 1980; 70: 419-439.
2. U.S. Department of Labor, Occupational Safety and Health Administration, 29 CFR Part 1910.1030. Occupational Exposure of Bloodborne Pathogenes; Final Rule. *Federal Register*; 56(235):64175, (1991).
3. USA Center for Disease Control/National Institute of Health Manual, Biosafety in Microbiological and Biomedical Laboratories", (1984).
4. National Committee for Clinical Laboratory Standards. Protection of Laboratory Workers from Instrument Biohazards and Infectious Disease Transmitted by Blood, Body Fluids, and Tissue: Approved Guideline. NCCLS Document M29-A, (1997).
5. Uotila, M., Ruoslahti, E. and Engvall, E., Two-site sandwich enzyme immunoassay with monoclonal antibodies to human alpha-fetoprotein. *J. Immunol. Methods*, 1981; 42:11-15).
6. Brizot, M.L., Jauniaux, E., Mckie, A.T., Farzaneh, F., and Nicolaides, K.H., Placental expression of α and β subunits of human chorionic gonadotrophin in early pregnancies with Down's syndrome. *Hum. Reprod.* 1995; 10: 2506-2509.
7. Forest, J., Masse, J., Rousseau, F., Moutquin, J., Brideau, N., and Belanger, M., Screening for Down Syndrome during the first and second trimesters: Impact of risk estimation parameters. *Clin. Biochem.* 1995; 28: 443-449.
8. Breimer, L., First trimester biochemical screening for trisomy 21: the role of free β -hCG and pregnancy associated plasma protein. *Ann. Clin. Biochem.*, 1995; 32: 233.
9. Loncar, K., Barnabei, V.M., and Larsen, J.W. Jr., Advent of maternal serum for Down syndrome screening. *Obstet. Gynecol. Surv.*, 1995; 50: 316-320.
10. Densem, J., and Wald, N.J., The stability of blood samples for the measurement of the free β subunit of chorionic gonadotrophin. *Prenat. Diagn.*, 1995; 15: 94-95.
11. Ozturk, M., Berkowitz, R., Goldstein, D., Bellet, D., Wands, J.R., Differential production of human chorionic gonadotropin and free subunits in gestational trophoblastic disease. *Am. J. Obstet. Gynecol.*, 1988; 158:193-198.
12. Wald, N.J., Cuckle, H.S., Densem, J.W., et al., Maternal serum screening for Down's syndrome in early pregnancy. *Br. Med. J.*, 1998; 297:883-887.
13. Hay, D.L., Placental histology and the production of human choriogonadotrophin and its subunits in pregnancy. *Br. J. Obstet. Gynaecol.*, 1988; 95: 1268-1275.
14. Macri, J.N., Kasturi, R.V., Krantz, D.A., et al., Maternal serum Down syndrome screening: Free β -protein is a more effective marker than human chorionic gonadotropin. *Am. J. Obstet. Gynecol.*, 1990; 163:1248-1253.
15. Ozturk, M., Bellet, D., Manil, L., et. al., Physiological studies of human chorionic gonadotropin (hCG), α hCG, and β hCG as measured by specific monoclonal immunoradiometric assays. *Endocrinology*, 1987;

120: 549-558.

16. Cole, L.A., Hartle, R.J., Laferla, J.J., et al., Detection of the free beta subunit of human chorionic gonadotropin (hCG) in cultures of normal and malignant trophoblast cells, pregnancy sera, and sera of patients with choriocarcinoma. *Endocrinology*, 1983; 113:1176-1178.
17. Gaspard, U.J., Reuter, A.M., Deville, J-L, et al., Serum concentration of human chorionic gonadotropin and its alpha and beta subunits. *Clin. Endocrinol. (OXF)*, 1980; 13: 319-329.
18. Tietz, N.W. ed., *Clinical Guide to Laboratory Tests*, 3rd Edition, W.B. Saunders, Co., Philadelphia, 1995: 134-135.

Plate Layout

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H