



EAGLE
BIOSCIENCES

Salivary Cortisol Ultrasensitive ELISA

Catalog Number:

COR32-K01 (1 x 96 Wells)

For Research Use Only. Not for use in diagnostic procedures

v. 16 (31 DEC 24)

EAGLE BIOSCIENCES, INC.

20A NW Blvd., Suite 112, Nashua, NH 03063
Phone: 617-419-2019 Fax: 617-419-1110

WWW.EAGLEBIO.COM



INTENDED USE

The Eagle Biosciences Salivary Cortisol ELISA Assay kit is designed and validated for the quantitative measurement of salivary Cortisol by ELISA. This kit features one 60 minute incubation for maximum assay sensitivity and precision as well as a saliva-like matrix assay buffer system containing a neutralizing agent for acidic or basic saliva samples.

For further information about this kit, its application or the procedures in this insert, please contact the Technical Service Team at Eagle Biosciences, Inc at www.EagleBio.com or at 866-411-8023.

ASSAY BACKGROUND

In 1966, Katz and Shannon using the Porter-Silber method were able to determine corticosteroid concentrations in saliva and were able to show that concentrations of corticosteroids in saliva, were related to blood concentrations. The advent of immunoassay made it possible to measure minute amounts of steroid hormones in blood. Subsequent modification of those assays allowed their measurement in saliva as well. These early assays, however, lacked validity due to matrix differences between serum and saliva, poor sensitivity, and cumbersome extraction methods. Recently, several papers have been published on the determination of salivary Cortisol under varying physiological conditions using more specific and sensitive EIA and ELISA methods.

Cortisol (hydrocortisone, compound F) is the principal glucocorticoid secreted by the adrenal cortex. Adrenal secretion of cortisol is modulated by a complex negative feedback mechanism involving the central nervous system, hypothalamus, pituitary and adrenals. ACTH released from the pituitary augments adrenal secretion of cortisol. In turn, increased levels of cortisol suppress pituitary secretion of ACTH while falling levels of cortisol are associated with rising levels of ACTH. Normally there is diurnal variation of cortisol with highest values measurable in the morning samples and lowest values obtained in the late afternoon. Cortisol levels rise independently of this circadian rhythm in response to stress or depression. Increased cortisol production is associated with Cushing's Syndrome and adrenal tumors while decreased production of cortisol is associated with adrenal insufficiency (Addison's disease) and adrenocorticotrophic hormone (ACTH) deficiency.

In blood 90% of the circulating cortisol is firmly bound to cortisol binding globulin (CBG), 7% is weakly bound to albumin and only 1-3% is free or unbound. In saliva the majority of cortisol occurs in the free or unbound form and enters the saliva via intracellular mechanisms. Numerous studies consistently report a high correlation between serum and saliva cortisol indicating that salivary cortisol levels clinically confirm levels of cortisol in serum.

ASSAY PRINCIPLES

This ultrasensitive cortisol saliva ELSIA is based on the competition principal and microplate separation. Cortisol calibrators of known concentration, unknown amounts of cortisol in saliva samples and a fixed amount of cortisol (analog) conjugated to horse-radish peroxidase (cortisol-HRP) compete for the binding sites with a rabbit-polyclonal-antiserum bound to GARGG (goat anti-rabbit gamma globulin) coated wells of a microplate. After incubation, unbound components are washed away, enzyme substrate solution is added, and a blue color is formed. This reaction is stopped with an acid solution to produce a yellow color. The optical density is then read at 450 nm. The amount of cortisol-HRP detected is inversely proportional to the amount of cortisol in a sample.



REAGENTS PROVIDED AND REAGENT PREPARATION

1. **GARGG Plate**- 1, 96-well plate (12x8 breakable strips) coated with goat anti-rabbit gamma globulin placed in a resealable foil bag. Sufficient for 78 singlicate or 39 duplicate sample measurements
2. **CALIBRATORS** – 7 vials, ready-to-use vials
 - CAL 0:** 0 ng/mL.....5.0 mL vial
 - CAL 1:** 0.1 ng/mL.....1.0 mL vial
 - CAL 2:** 0.3 ng/mL.....1.0 mL vial
 - CAL 3:** 1.0 ng/mL.....1.0 mL vial
 - CAL 4:** 3.0 ng/mL.....1.0 mL vial
 - CAL 5:** 10.0 ng/mL.....1.0 mL vial
 - CAL 6:** 30.0 ng/mL.....1.0 mL vial
3. **CONTROL 1 (LOW)** – 1 vial of 1.0 mL, concentration on vial, Range on QC
4. **CONTROL 2 (HIGH)** – 1 vial of 1.0 mL, concentration on vial, Range on QC
5. **ANTI-CORTISOL ANTIBODY** – 1 vial of 6 ml of antibody produced in rabbit. Diluted in phosphate buffer base. Contains animal protein and a binding protein blocker in blue solution.
6. **HRP CONCENTRATE (10X)** – 1 amber glass vial of 0.7 mL
7. **HRP CONJ BUFFER** – 6.3 mL of yellow solution to be used for working reagent preparation only
8. **HRP WORKING REAGENT** – Preparation;
Determine the amount of working cortisol-HRP needed and dilute 1:10 in conjugate buffer (# 7). For example, mix 0.5 mL of cortisol-HRP concentrate (#6) + 4.5 mL of Cortisol-HRP conjugate buffer (#7). This is sufficient for 100 wells. Immediately after use, store the unused portion of the Cortisol-HRP working solution at 2-8°C. Discard if not used within 4 (four) weeks of mixing.
9. **WASH SOLUTION (10x)** – 1 vial of 50 mL. Dilute 1:10 with DI Water
10. **COLOR DEVELOPMENT REAGENT (TMB)** – 1 vial of 12 mL of TMB plus H₂O₂. Light sensitive
11. **STOP SOLUTION** – 1 vial of 12 mL acid solution

*Concentration of cortisol calibrators and controls are actual and traceable to NIST Cortisol Catalog #SRM921 Lot#921

STORAGE AND STABILITY

- When stored at 2-8°C, unopened reagents will retain activity until the expiration date. Do not use reagents beyond this date
- Use only reagents supplied with this kit. Do not interchange reagents with different lot numbers
- Opened reagents must be stored at 2-8°C
- Microtiter wells must be stored at 2-8°C. Once foil bag has been opened, care should be taken to reseal tightly.
- Opened kits retain activity one (1) month if stored properly
- Expiration dates and lot numbers are printed on the labels

MATERIALS NEEDED BUT NOT SUPPLIED

1. Device to dispense very accurately 25 µl of saliva.
2. Multichannel pipettors.
3. Microplate or orbital shaker
4. Vortex Mixer
5. Microplate washer (not required, plates can be washed manually).
6. Microplate reader capable of reading 450 nm with 4 parameter data reduction or comparable software.
7. Plate sealer
8. Collection device



SAMPLE COLLECTION AND PREPERATION

Sample collection and processing procedure must be followed.

1. Avoid food consumption, drinking coffee or alcohol, smoking, or chewing gum one (1) hour prior to sample collection.
2. Rinse mouth thoroughly with water 15 minutes prior to collection
3. Collect whole saliva by unstimulated passive drool by allowing saliva to drip off the lower lip into a graduated plastic tube or by allowing saliva to accumulate in the floor of the mouth and spitting it into the recommended collection device.
4. Time and date specimen. Refrigerate then freeze (-20C or below) samples until day of assay. On day of assay, thaw samples to facilitate precipitation of mucins. Centrifuge at 1500xg for ten (10) minutes. Bring samples to room temperature and assay

Sample Stability

STORAGE	Room Temperature 20-30°C	37°C	2-8°C	≤ -15°C (7 freeze / thaw cycles)	≤ -15°C (Long term)
STABILITY	Up to 7 days	Up to 7 Days	Up to 7 Days	Up to 7 Days	Up to 180 Days

Interferences

An in-vitro experiment was performed by spiking three (3) levels of cortisol with high concentrations (> 1000 fold those of cortisol) of five (5) commonly consumed products; alcohol, coffee (as caffeine), cigarette (as nicotine) and food and gum as extracts. The results obtained demonstrate no significant difference between the controls and the spiked samples.

Calibrator Control Sample (I.D.)	Calibrator control (µL)	Cortisol-HRP Working Reagent (µL)	Anti-Cortisol (µL)		Diluted 10X Wash Solution (µL)		Color Developer (µL)		Stopping Solution (µL)	
0	25	50	50	Mix. Incubate 60 min. at room temperature. with shaking	300	Wash 3x	100	Mix. Incubate 30 min. At room temperature	100	Mix. Read at 450nm
0.1	25	50	50		300		100			
0.3	25	50	50		300		100			
1	25	50	50		300	100				
3	25	50	50		300	100				
10	25	50	50		300	100				
30	25	50	50		300	100				
Control #1	25	50	50		300	100				
Control #2	25	50	50		300	100				
Sample #1	25	50	50		300	100				



ASSAY PROCEDURE

1. To GARGG microplate pipet **25 µL** of ready-to-use Salivary Cortisol ELISA calibrators (0, 0.1, 0.3, 1.0, 3.0, 10.0 and 30 ng/mL), controls, and saliva samples.
2. Add **50 µL** of Cortisol-HRP working reagent to all wells (see Reagent Preparation Section-number 8).
3. Add **50 µL** of Cortisol ELISA antibody.
4. Cover microplate with plastic sealer. Incubate by shaking on a microplate orbital shaker set at 500-900 rpm for **60 minutes** at room temperature.
5. After incubation, decant the contents of the wells. **Wash** 3 times with **300 µL** of diluted wash solution (10 mL of 10X Wash solution ELISA #1 diluted with 90 mL of D.I. water). After the 3rd wash, invert GARGG microplate on an absorbent paper and tap dry.
6. Pipette **100 µL** of TMB into each well. Shake briefly (manual). **Incubate** for 30 **minutes** at room temperature.
7. Pipette **100 µL** of Stopping Solution ELISA #1 into each microtiter well of the GARGG plate. Shake briefly (manual). Color changes from blue to yellow.
8. **Read** at 450 nm on a microplate reader within **30 minutes**.

Note: If samples exceed the highest calibrator, dilute with zero calibrator and make appropriate concentration correction.

TYPICAL RESULTS

Typical Calibration Curve for 25 µL Sample Size			
Calibrators (ng/L)	Mean Absorbance (450 nm)	%B/B0	Value (ng/mL)
0	2.231	100.0	0.0
0.1	2.056	92.2	0.1
0.3	1.717	77.0	0.3
1	1.155	51.8	1.0
3	0.614	27.5	3.0
10	0.246	11.0	10
30	0.130	5.8	30
Control #1(low)	1.130	50.6	1.04
Control #2 (high)	0.159	7.1	20.14
Sample #1	0.970	43.5	1.40
Sample #2	0.235	10.5	11.26

Conversion factor ng/mL to nmol/L=2.76

CALCULATION

1. Compute the average optical density (OD) for the zero (B0) calibrator
2. Calculate the percent bound (B/B0) for each calibrator, control and unknown by dividing the average OD (B) by the average OD for the zero (B0) x 100
3. Plot percent bound (B/B0) versus the calibrator concentrations and draw the best fit for the curve.
4. Plot percent bound (B/B0) of the controls and unknowns to determine saliva cortisol concentrations.
5. Alternatively, determine the concentrations of the controls and unknowns by interpolation using software capable of logistics using a 4-parameter sigmoid minus curve fit.

Analytical measuring range (AMR): 0.1 ng/ml-30.0 ng/ml



Samples with cortisol values greater than 30 ng/ml (82.77 nmol/L) should be diluted 1:10 with zero (0) calibrator and rerun for accuracy. Obtain the final cortisol concentration by multiplying the diluted sample by the dilution factor.

QUALITY CONTROL

The expected values for the controls are stated on the label of each control which are included in the kit. The results can only be accepted if the expected values are met.

EXPECTED VALUES

AM Expected Values				
Subjects (Number)	Subjects (Gender)	Age (Years)	AM Median (ng/mL)	AM Range (ng/mL)
152	76 Males 76 Females	23-68	6.70	2.58-12.69

PM Expected Values				
Subjects (Number)	Subjects (Gender)	Age (Years)	AM Median (ng/mL)	AM Range (ng/mL)
152	76 Males 76 Females	23-68	0.58	0.25-2.96

*it is recommended that each laboratory establishes its own range of normal values.

PERFORMANCE CHARACTERISTICS

Specificity of Antiserum

Compounds	Spiked Concentration (ng/mL)	% Cross-reactivity
C-21 Steroids		
Cortisol	10,000	100.0
17-OH-Progesterone	10,000	0.0284
Pregnenolone	10,000	0.0038
17-OH-Pregnenolone	10,000	0.0066
Progesterone	10,000	0.0079
Deoxycorticosterone	10,000	0.0517
11-Desoxycortisol	10,000	1.8133
Dexamethasone	10,000	0.0164
Cortisone	10,000	0.7600
Corticosterone	10,000	1.0847
Aldosterone	10,000	0.0070
C-19 Steroids		
Androstenedione	10,000	0.0038
Testosterone	10,000	0.0042
5αDHT	10,000	0.0019
DHES-SO ₄	10,000	0.0031
Androstanedione	10,000	0.0028
C-18 Steroids		
Estradiol 17β	10,000	0.0024

Estradiol 17a	10,000	0.0003
Estrone	10,000	0.0010
Estriol	10,000	0.0015
Other structurally related steroids		
Dehydroisoandrosterone	1000	0.0076
6methyl-17-hydroxyprogesterone	1000	0.1427
6β-Hydroxycortisol	1000	1.7177
Prednisone	1000	1.0874
Prednisolone	1000	25.9001

At >10% cross reaction prednisolone is a potential substance

Detection Limit

The LOB (limit of the blank), the LOD (limit of detection) and the LOQ (limit of quantitation) were determined by generating one hundred twenty (120) measurements of cortisol free saliva and low level (<0.1 ng/mL) cortisol samples

Limit of the Blank (LOB) ng/mL	Limit of Detection (LOD) ng/mL	Limit of Quantitation (LOQ) ng/mL
0.0392	0.0519	0.0519

Precision and Reproducibility

Intra-assay

The intra-assay precision was determined from the mean of 20 replicates of low, medium, and high samples

Sample	N	Mean (ng/mL)	Standard Deviation (ng/mL)	%CV
Low	20	0.627	0.034	5.4
Medium	20	3.995	0.266	6.7
High	20	25.232	1.579	6.3

Inter-assay

The inter-assay precision was determined from the mean of the average duplicates of 12 separate assays with low, medium and high samples

Sample	N	Mean (ng/mL)	Standard Deviation (ng/mL)	%CV
Low	12	0.587	0.037	6.3
Medium	12	4.163	0.301	7.2
High	12	25.126	0.712	2.8

Repeatability

This study was conducted during 4 days of a familiarization period and 20 days of testing. Two assays were performed daily with a minimum of 2 hours between assays, three (3) different reagent lots and three (3) saliva pools were used for the study (low, medium, and high). The pools were aliquoted and frozen until day of assays.

Precision Low Concentration Pool

	Standard Deviation, (SD)	% Coefficient of Variation (CV)
Within Run	0.0224	3.79
Between Run	0.0426	7.80
Repeatability	0.0162	2.73
Total Device Precision	0.0538	9.09

Precision Medium Concentration Pool

	Standard Deviation, (SD)	% Coefficient of Variation (CV)
Within Run	0.1475	3.60
Between Run	0.0514	1.26
Repeatability	0.1025	2.50
Total Device Precision	0.1869	4.56

Precision High Concentration Pool

	Standard Deviation, (SD)	% Coefficient of Variation (CV)
Within Run	0.4442	1.76
Between Run	0.2915	1.15
Repeatability	0.62276	2.46
Total Device Precision	0.8185	3.24

Inter-lot Variation

The inter-lot precision was determined by duplicate measurements of five (5) saliva pools and three (3) spiked controls in saliva like matrix.

Saliva Samples	Lot #012	Lot #013	Lot #014	Inter-lot	Inter-lot	Inter-lot
ID	Mean (ng/ml)	Mean (ng/ml)	Mean (ng/ml)	Mean (ng/ml)	SD (ng/ml)	CV (%)
20	4.65	4.45	4.79	4.64	0.164	3.5
21	0.67	0.61	0.71	0.67	0.049	7.4
22	2.02	1.95	2.09	2.02	0.069	3.4
23	4.75	4.69	4.76	4.73	0.041	0.9
24	2.01	1.99	2.04	2.01	0.026	1.3
25	3.64	3.67	3.71	3.68	0.036	1
LC	0.98	0.94	1.01	0.98	0.036	3.7
MC	5.21	5.31	5.49	5.34	0.14	2.6
HC	10.79	10.13	10.52	10.48	0.329	3.1

Linearity Study

Ten (10) sample concentrations that span the assay measuring range were performed per EP6-A, Evaluation of the Linearity of Quantitative Measurement Procedures

S= 10 samples (dilutions)

$$\text{Concentration} = (C1 \cdot V1 + C10 \cdot V10) / (V1 + V10)$$



	C1 (ng/mL)	V1 (ng/mL)	C10 (ng/mL)	V10 (ng/mL)	Calculated Concentration (ng/mL)	Obtained Concentration (ng/mL)	Recovery (%)
1	0.093				0.100	0.093	93
2	0.093	0.889	33.788	0.111	3.833	3.729	97.3
3	0.093	0.778	33.788	0.222	7.573	7.620	100.6
4	0.093	0.667	33.788	0.333	11.313	10.842	95.8
5	0.093	0.556	33.788	0.444	15.054	14.350	95.3
6	0.093	0.444	33.788	0.556	18.827	18.313	97.3
7	0.093	0.333	33.788	0.667	22.568	21.547	95.5
8	0.093	0.222	33.788	0.778	26.308	24.694	93.9
9	0.093	0.111	33.788	0.889	30.048	30.459	101.4
10					35.000	33.788	96.6

Recovery

Ten (10) saliva samples containing different levels of endogenous cortisol were spiked with known quantities of cortisol and assayed.

Sample	Endogenous (ng/mL)	Added (ng/mL)	Expected (ng/mL)	Observed (ng/mL)	Recovery (%)
1	0.493	0.250	0.743	0.739	99.5
2	0.878	0.500	1.378	1.291	93.7
3	1.551	1.000	2.551	2.641	103.5
4	1.850	2.000	3.850	3.958	102.8
5	0.936	4.000	4.936	4.951	100.3
6	1.042	8.000	9.042	9.394	103.9
7	0.691	16.000	16.691	17.165	102.8
8	0.622	20.000	20.622	19.997	97.0
9	2.057	24.000	26.057	24.938	95.7
10	0.348	28.000	28.348	28.943	102.1

Method Comparison

A comparative study was performed between this assay and an FDA cleared predicted device. A total of 160 samples were used for the study of which 6 samples were spiked representing 3.75% of the total number of samples. The results show the following regression and correlation statistics

Linear Regression Equation: $Y=1.0269X + 0.0994$

Correlation: $R^2=0.989$

LIMITATIONS

- The reagents are optimized to measure cortisol directly in saliva
- Cortisol levels are elevated during the later stages of pregnancy and in women on contraceptives or after long-term use of contraceptives (28, 29)
- Elevated cortisol levels can be found in conditions of sepsis, infection, chronic liver disease, and renal failure. Low cortisol levels result from liver disease, pituitary hyposecretion, hypothyroidism, or steroid therapy
- Note that an in-vitro study to identify potential interfering substances in the measurement of salivary cortisol, may not identify some interferents and the form(s) of potential interferents being tested may not represent the naturally occurring forms.



- The use of topical creams or ointments containing hydrocortisone (true cortisol) should be avoided as they can cause preanalytical contamination of the saliva sample indistinguishable from endogenous cortisol as measured by immunoassay or LC-MS/MS (30).

PRECAUTIONS

- This kit is for research use only
- Compare contents and packing list, if there is breakage or shortage, notify Eagle Biosciences immediately
- Do not pipette reagents by mouth
- Do not smoke, eat or drink while performing assay
- Wear disposable gloves and proper lab protection and attire
- Treat all samples as potentially infectious
- Do not mix reagents from other lots
- Avoid contact with TMB and Stop solutions. If contact occurs, wash thoroughly with water
- Eagle Biosciences is not responsible for outcomes as results of tampering with the reagents or using them unconventionally

REFERENCES

1. Katz FH, Shannon IL: Parotid fluid corticosteroids a good measure of cortisol available to the cell: *J.Clin.Invest.* 45:1031, 1966.
2. Katsh FH, Shannon IL: Parotid fluid cortisol and cortisone: *J.Clin.Invest.* 48:448, 1969.
3. Katz FH, Shannon IL: Adrenal corticosteroids in submaxillary fluid: *J. Dent. Res.* 48:448, 1969.
4. Filare E, Duche P, Lac G, Robert A: Saliva cortisol, physical exercise and training: influences of swimming and handball on cortisol concentrations in women: *Eur. J. Appl. Physiol.* 74: 274, 1996.
5. Filare E, Duche P, Lac G: Effect of training for two ball games on the saliva response of adrenocortical hormones to exercise in elite sports women: *Eur.J.Appl.Physiol.* 77:452, 1998.
6. Filare E, Duche P, Lac G: Effects of amount of training on the saliva concentrations of cortisol, Dehydroepiandrosterone: Cortisol concentration ration in women over 16 weeks of training. *Euro. J. Appl. Physiol.* 78:466, 1998.
7. Aardal-Erikson E, Karlberg BE, Holm AC: Salivary cortisol-an alternative to serum cortisol determinations in dynamic function tests: *Clin. Chem. Lab, Med.*, 4:15, 1998.
8. Smyth J, Ochenfels MC, Porter L, Kirschbaum C, Hellhammer DH, Stone A, *Psychoneuroendocrinology* 23:353, 1998.
9. Chicharro JL, Lucia A, Perez M, Vaquero AF, Urena R: Saliva composition and exercise: *Sports Med.* 26:17, 1998.
10. Grassi G, Ahs A, Lundberg U: Psychological correlates of salivary cortisol secretion amount unemployed men and women: *Integr. Physiol. Behav. Sci.*: 23:249, 1998.
11. Ratt H, Ratt JL, Findling JW: Late-night salivary cortisol as a screening test for Cushing's Syndrome: *J. Clin. Endocrinol. Metab.* 83:2681, 1996.
12. Findling JW and Ratt H: newer diagnostic techniques and problems in Cushing's Disease: *endocrinol. Metab. Clin. North Am.* 28:19, 1999
13. Van Honk J, Tuiten A, Van der Hout M, Koppeschaar H, Thijssen J, de Hann E, Verbaten R: Baseline salivary cortisol levels and preconscious selective attention for threat a pilot study: *Psychoendo.* 23, 741, 1996.
14. Odler J, Cawood EH, Bancroft J: Salivary cortisol in women with and without perimenstrual mood changes; *J. Psychosom, Res.* 45:557, 1998.
15. Raff H, Ratt JL, Duthie Eh, Wilson CR, Sasse EA, Rudman I, Mattison D: Elevated Salivary Cortisol in the evening in health elderly men and women correlation with bone mineral density. *J. Gerontol A Biol. Sci. Med, Sci.* 54:M479, 1999.
16. Aardal-Eriksson E, Eriksson TE, Holm AC, Lundin T,: Salivary cortisol and serum prolactin in relation to stress rating scales in a group of rescue workers. *Biol. Psychiatry* 45:850, 1999.



17. Filare E, Scatt C, Duche P, Lac G: The relationship between salivary adrenocortical hormones changes and personality in elite female athletes during handball and volleyball competition: *ResQ, Exerc. Sport*: 70:297, 1999.
18. Aardel E and Holm AC. Cortisol in saliva reference ranges and relation to cortisol to serum. *European J of Clinical Chemistry and Clinical Biochemistry* 33: 927, 1995.
19. Clements Ad and Parker CR: The relationship between salivary cortisol concentrations in frozen versus mailed samples. *Psychoendocrinology* 23:613, 1998.
20. Schwartz EB, Granger DA, Susman EJ, Gunnar MR and Laird B: Assessing salivary cortisol in studies of child development., *Child development* 69, 1503-1513, 1998.
21. Korenman, S.G.,D.K. Grammer and B.M. Sherman, *Practical Diagnosis: Endocrine Disease*. Houghton Mitten Professional Publishers, 1978.
22. Harshman, J.M., *Endocrine Pathophysiology: A Patient Oriented Approach*, Lea and Fabiger, 1977.
23. Garza, R.and G.E. Abraham, *Analytical Letters*, 5:767, 1972.
24. Abraham, G.E., Ed., *Radioassay Systems in Clinical Endocrinology* Marcel Dekker, Inc. New York, 1961.
25. Vining, R.F., McGinley, R.A. and Symons, R.G., (1983). Hormones in Saliva mode of Entry and Consequent Implications for Clinical Interpretation. *Clin. Chem.*, 29, 1752-1756.
26. Francis, S.J. et. al. (1987) Assesment of Andrenocortical activity in term newborn infants using salivary cortisol determinations. *J. of Pediatrics*, 111, 129-333.
27. Hiramtsu, R. (1981) Direct Assay of Cortisol in human saliva by solid phase radioimmunoassay and its clinical applications. *Clinica Chimica. ACTA.* 177, 239-240.
28. Vining, R.F., et al. (1983). Salivary Cortisol: A better measure of adrenal cortical function than serum cortisol. *Annals Clin. Biochem*, 20, 329-335.
29. Susman, E.J., et. al. (1999). Corticotropin-releasing hormone and cortisol: Longitudinal Associations with depression and antisocial behavior in pregnant adolescents. *J. of the American Academy of Child and Adolescent Psychiatry*, 38, 460-467.
30. Raff, H., and Ravinder, J.S. Measurement of Late-Night Salivary Cortisol and Cortisone by LC-MS/MS to assess preanalytical sample contamination with topical hydrocortisone. *Clin. Chem.* 2012; 58: 947-48

WARRANTY INFORMATION

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.