



REFERENCE RANGES

Reference ranges (95%) were estimated using samples obtained from individuals of diverse races. Each laboratory shall establish their own range of reference values.

Serum Cohort Group	n	Median (ng/mL)	95% Range (ng/mL)	Total Range (ng/mL)
Adult Males and Non-Pregnant Females	120	< 0.058	ND–0.11	ND–0.12
Pregnant Females First Trimester	30	0.15	—	ND–2.95
Pregnant Females Second Trimester	50	1.20	0.46–3.04	0.45–3.07
Pregnant Females Third Trimester	25	9.5	—	3.6–14.3

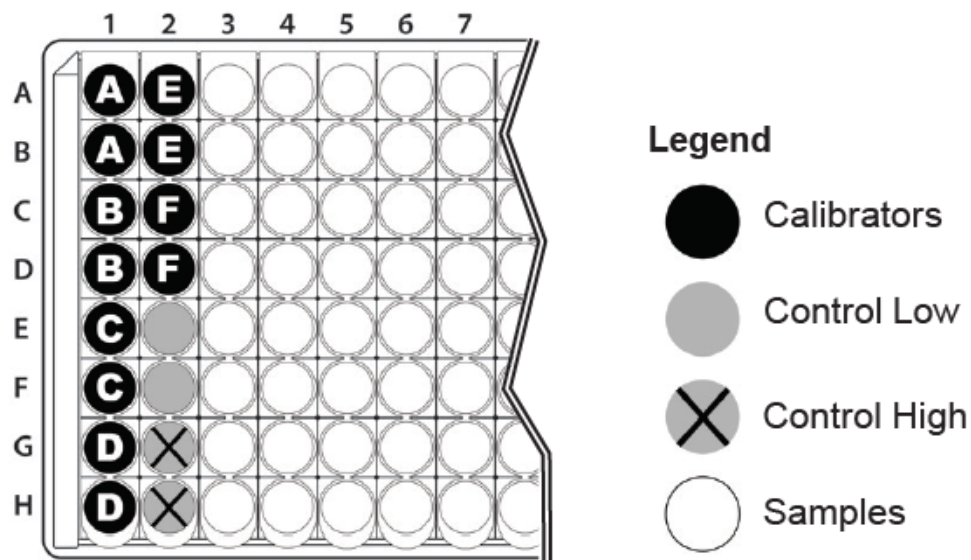
ND = Non-determined; value less than the limit of detection.



LITERATURE

1. Berkane N, et al. From Pregnancy to Preeclamsia: A Key Role for Estrogens. *Endocr Rev.* 2017; 38(2):123-144.
2. Fleck SC, et al. Comparative estrogenicity of endogenous, environmental and dietary estrogens in pregnant women I: Serum levels, variability and the basis for urinary biomonitoring of serum estrogenicity. *Food Chem Toxicol.* 2018;115:511-522.
3. Kos-Kudla B, et al. Comparative studies of serum and salivary estriol concentrations in third trimester of normal pregnancy. *Med Sci Monit.* 1999;5(2):285-288.
4. Falah N, et al. Estriol review: Clinical applications and potential biomedical importance. *Clin Res Trials.* 2015;1(2):29-33.
5. Johnsson VL, et al. Plasma progesterone, estradiol, and unconjugated estriol concentrations in twin pregnancies: Relation with cervical length and preterm delivery. *Acta Obstet Gynecol Scand.* 2019;98(1):86-94.
6. Montazeri S, et al. Investigating Relationship between Saliva Estriol to Progesterone Ratio and Preterm Delivery. *International Journal of Pharmaceutical Research & Allied Sciences.* 2016;5(3):264-9.
7. Hampson E, et al. Steroid concentrations in antepartum and postpartum saliva: normative values in women and correlations with serum. *Biol Sex Differ.* 2013;4(1):7.

RECOMMENDED MICROPLATE LAYOUT





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.