

## Progesterone Metabolites Competitive ELISA Kit

Catalog Number EIAPROGM (96 tests)

Pub. No. MAN0018783 Rev A.0

**Note:** For safety and biohazard guidelines, see the “Safety” appendix in the *ELISA Technical Guide* (Pub. no. MAN0006706). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

### Product description

The Progesterone Metabolites ELISA Kit is designed to detect and quantify progesterone metabolites in dried fecal extracts and urine samples. The assay recognizes progesterone metabolites independent of species.

Progesterone is a C-21 steroid hormone that is involved in the female menstrual cycle, gestation, and embryogenesis of humans and other species. In different animal species, progesterone can be metabolized and excreted as a variety of progesterone molecules. Measurement of these progesterone molecules can provide vital data to aid reproductive strategies in endangered species.

### Contents and storage

Kit and components are shipped at -20°C. Upon receipt, store the kit at -20°C. Once open, store the kit at 4°C and use within 2 weeks.

Components	Quantity
Coated Clear 96-Well Plate, clear plastic microtiter plate coated with goat anti-rabbit IgG	1 plate
Progesterone Standard; 1,000 ng/mL	40 µL
Progesterone Antibody	3 mL
Progesterone Conjugate	3 mL
Assay Buffer Concentrate (5X)	28 mL
Wash Buffer Concentrate (20X)	30 mL
TMB (Tetramethylbenzidine) Substrate	11 mL
Stop Solution; contains 1 M HCl, <b>CAUSTIC</b>	5 mL
Plate Sealer	1

### Materials required but not supplied

- Distilled or deionized water
- Microtiter plate reader with software capable of measurement at or near 450 nm (preferably with correction between 570 nm and 590 nm).
- Microplate shaker.
- Plate washer—automated or manual (squirt bottle, manifold dispenser, or equivalent)
- Calibrated adjustable precision pipettes and glass or plastic tubes for diluting solution
- Ethanol for extraction of fecal samples.

### Prepare 1X Wash Buffer

1. Dilute Wash Buffer Concentrate 1:20 by adding 1 part of the Concentrate to 19 parts of deionized or distilled water. Label as 1X Wash Buffer.
2. 1X Wash Buffer is stable at room temperature for 3 months.

### Prepare 1X Assay Buffer

1. Dilute Assay Buffer Concentrate 1:5 by adding 1 part of the Concentrate to 4 parts of deionized or distilled water. Label as 1X Assay Buffer.
2. Store the Concentrate and 1X Assay Buffer in the refrigerator. 1X Assay Buffer is stable at 4°C for 3 months.

### Procedural guidelines

**IMPORTANT!** Reagents are lot-specific. Do not mix or interchange different reagent lots from various kit lots.

- Review the **Procedural guidelines** and **Plate washing directions** in the *ELISA Technical Guide* available at [thermofisher.com](http://thermofisher.com).
- Allow reagents to reach room temperature before use. Mix to redissolve any precipitated salts.
- Solutions containing sodium azide will inhibit the activity of the peroxidase conjugate. Ensure that there is no contamination of labware or the plate washer with azide-containing solutions.

**For research use only. Not for use in diagnostic procedures.**

## Sample preparation guidelines

- Refer to the *ELISA Technical Guide* at [thermofisher.com](http://thermofisher.com) for detailed sample preparation procedures.
- Collect samples in pyrogen/endotoxin-free tubes.
- Freeze samples after collection if samples will not be tested immediately. Avoid multiple freeze-thaw cycles of frozen samples. Thaw completely and mix well (do not vortex) prior to analysis.
- If large amounts of particulate matter are present in the sample, centrifuge or filter sample prior to analysis.

## Prepare samples

Sample concentrations should be within the range of the standard curve. Because conditions may vary, each investigator should determine the optimal dilution for each application.

Use all samples within **2 hours** of dilution, or store at  $-20^{\circ}\text{C}$  or lower until ready to perform assay.

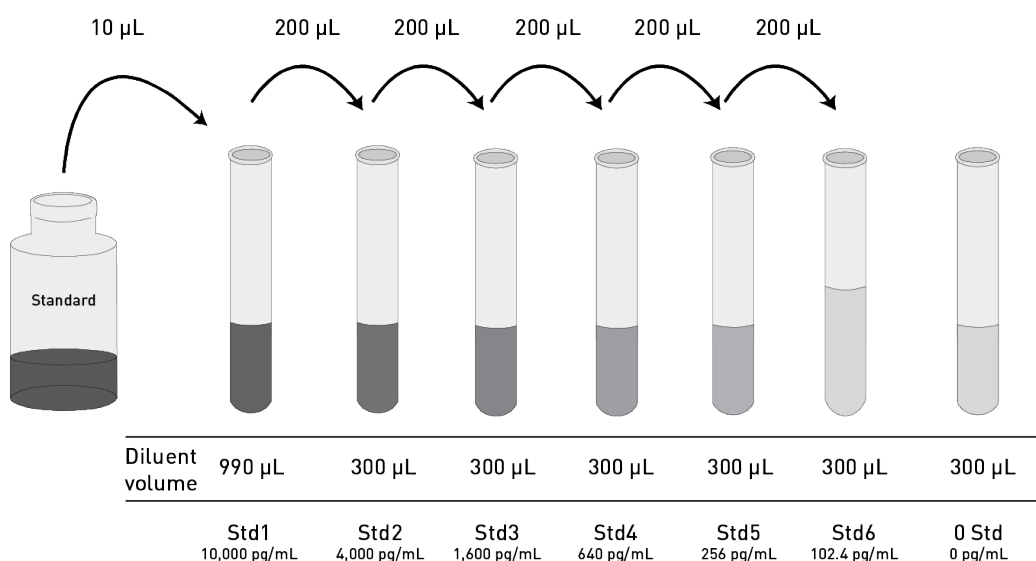
Sample type	Procedure
Urine	Dilute samples $\geq 1:8$ with 1X Assay Buffer. <b>Note:</b> A Urinary Creatinine Detection Kit (Cat. No. EIACUN) is available for measuring urine creatinine for normalization of progesterone metabolites in a random urine specimen.
Dried feces	See detailed extraction protocol on the product page at <a href="http://thermofisher.com">thermofisher.com</a> <b>Note:</b> The ethanol concentration in the final diluted Assay Buffer dilution added to the well should be $<5\%$ .

## Dilute standards

**Note:** Use glass or plastic tubes for diluting standards.

The Progesterone Standard contains an organic solvent. Pipette the standard up and down several times to wet the pipet tip before transfer to ensure that volumes are accurate.

1. Add 10  $\mu\text{L}$  Progesterone Standard to one tube containing 990  $\mu\text{L}$  1X Assay Buffer and label as 10,000 pg/mL Progesterone.
2. Add 300  $\mu\text{L}$  1X Assay Buffer to each of 6 tubes labeled as follows: 4,000, 1,600, 640, 256, 102.4, and 0 pg/mL Progesterone.
3. Make serial dilutions of the standard as described below in the dilution diagram. Mix thoroughly between steps.
4. **Use the standards within 2 hours of preparation.**



## Perform ELISA (Total assay time: 1.5 hours)

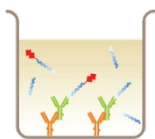
**IMPORTANT!** Perform a standard curve with each assay.

Allow all components to reach room temperature before use. Mix all liquid reagents prior to use.

Determine the number of 8-well strips required for the assay. Insert the strips in the frames for use. Re-bag any unused strips and frames, and store desiccated at 2°C to 8°C for future use. The silica pack in the bag keeps the plate dry, and turns from blue to pink if the bag is not properly sealed.

### Bind antigen

- Add 50 µL of standards or samples (see "Prepare samples" on page 2) to the appropriate wells.
- Add 75 µL of 1X Assay Buffer to wells for detecting non-specific binding (NSB).
- Add 50 µL of 1X Assay Buffer to wells for detecting maximum binding (B0 or zero standard)
- Add 25 µL of Progesterone Conjugate to each well.
- Add 25 µL of Progesterone Antibody to each well except NSB wells.
- Tap the side of the plate to mix. Cover the plate with plate sealer.
- Shake at room temperature for 1 hour. If the plate is not shaken signals bound will be approximately 25% lower.
- Thoroughly aspirate the solution and wash wells 4 times with 300 µL of 1X Wash Buffer.



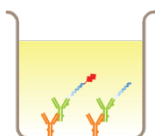
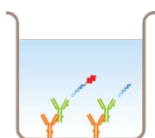
### Add chromogen

- Add 100 µL TMB Substrate to each well. The substrate solution will begin to turn blue.
- Incubate for 30 minutes at room temperature without shaking.

**Note:** TMB should not touch aluminum foil or other metals.

### Add stop solution

Add 50 µL Stop Solution to each well. Tap side of the plate gently to mix. The solution in the wells changes from blue to yellow.



## Read the plate and generate the standard curve

- Read the absorbance at 450 nm. Read the plate within 10 minutes after adding the Stop Solution.
- Average the duplicate Optical Density (OD) values for each standard and sample. Use curve-fitting software to generate the standard curve. A four parameter algorithm provides the best standard curve fit. Optimally, the background absorbance may be subtracted from all data points, including standards, unknowns and controls, prior to plotting.
- Calculate the concentrations for unknown samples and controls from the %B/B0 curve. Multiply value(s) obtained for sample(s) by the appropriate factor to correct for the sample dilution.

**Note:** Dilute samples producing signals lower than that of the highest standard in 1X Assay Buffer and reanalyze. Multiply the concentration by the appropriate dilution factor.

## Performance characteristics

### Standard curve (example)

The following data were obtained for the various standards over the range of 0–10,000 pg/mL Progesterone.

Progesterone Standard (pg/mL)	Net Optical Density (450 nm)*	%B/B0
10,000	0.132	10.3
4,000	0.258	20.2
1,600	0.472	37.1
640	0.755	59.3
256	0.964	75.7
102.4	1.128	88.6
0	1.273	100

**Note:** The NSB gave a Mean OD result of 0.078.

### Intra-assay precision

Samples were assayed in replicates of 20 to determine precision within an assay.

Parameters	Sample 1	Sample 2	Sample 3
Mean (pg/mL)	2,377	1,517	710.3
%CV	4.9	7.2	8.0

CV = Coefficient of Variation

### Inter-assay precision

Samples were assayed in duplicates in 19 assay runs by six operators to determine precision between assays.

Parameters	Sample 1	Sample 2	Sample 3
Mean (pg/mL)	2,442	1,665	844.6
%CV	6.8	6.6	8.8

CV = Coefficient of Variation

## Performance characteristics, continued

### Expected values

Four human urine samples were tested in the assay, three of which came from pregnant women who were 17, 21, and 26 weeks pregnant. Adjusted neat concentrations of progesterone ranged from 39.46 to 4312.6 ng/mL. When adjusted for urine creatinine using the Urinary Creatinine Detection Kit (Cat. No. EIACUN), the values ranged from 49.7 to 215.9 ng/mg creatinine.

### Sensitivity

The minimum detectable dose of Progesterone is 51.2 pg/mL. This was determined by adding two standard deviations to the mean O.D. obtained when the zero standard was assayed 20 times, and calculating the corresponding concentration.

### Linearity

Linearity was determined for diluted urine by taking two samples, one with a low level (179.2 pg/mL) and one with a higher level (1,729 pg/mL) of epiandrosterone, and mixing them in the ratios given below. The measured concentrations were compared to the expected values based on the ratios used.

#### Urine

Low Sample %	High Sample %	Expected Conc. (pg/mL)	Observed Conc. (pg/mL)	% Recovery
80	20	489.1	519.9	106.3
60	40	799.1	797.5	99.8
40	60	1,109	1,174	105.9
20	80	1,419	1,451	102.2

Mean Recovery 103.6%

### Specificity

The following cross reactants were tested in the assay and calculated at the 50% binding point.

Steroid	Cross-reactivity %
Progesterone	100.00
5 $\beta$ -dihydroprogesterone	61.90
5 $\alpha$ -dihydroprogesterone	56.70
Pregnanolone (5 $\beta$ -pregnan-3 $\alpha$ -ol-20-one)	41.20
Epiallopregnanolone (5 $\alpha$ -pregnan-3 $\beta$ -ol-20-one)	38.30
Allopregnanolone	27.30
Pregnenolone	17.60
Epipregnanolone	10.20
17 $\alpha$ -hydroxyprogesterone	5.70
11 $\alpha$ -hydroxyprogesterone	4.90
20 $\alpha$ -hydroxyprogesterone	0.34
Allopregnandiol	0.29

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



**Manufacturer:** Life Technologies Corporation | 7335 Executive Way | Frederick, MD 21704 | USA

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

© 2019 Thermo Fisher Scientific Inc. All rights reserved. MAGPIX, FLEXMAP 3D, xPONENT, Luminex, Luminex 100, and Luminex 200 are trademarks of Luminex Corporation. Excel is a trademark of Microsoft Corporation. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.