



**EAGLE**  
BIOSCIENCES

# Melatonin Saliva ELISA Assay Kit

Catalog Number: MEL32-K01

For Research Use Only. Not for use in diagnostic procedures.  
v. 6 (31 DEC 22)

---

EAGLE BIOSCIENCES, INC.  
20A Northwest Blvd., Suite 112, Nashua, NH 03063  
Phone: 617-419-2019 Fax: 617-419-1110  
[WWW.EAGLEBIO.COM](http://WWW.EAGLEBIO.COM)



## INTENDED USE

The Eagle Biosciences Melatonin Saliva ELISA Assay Kit is designed and validated for the direct quantitative measurement of melatonin in human saliva. The Melatonin saliva ELISA Assay Kit is for research use only and not to be used for diagnostic procedures.

## ASSAY BACKGROUND

Melatonin (N-Acetyl-5-methoxytrptamine) is a biogenic amine that is found in animals and plants. In mammals, melatonin is produced by the pineal gland. Its secretion increases in darkness and decreases during exposure to light. Melatonin is implicated in the regulation of sleep, mood and reproduction. Melatonin is also an effective antioxidant.

## PRINCIPLE OF THE ASSAY

The Eagle Biosciences Direct Salivary Melatonin EIA kit is based on the competition principal and microplate separation. Melatonin in standards (calibrators) and samples compete with a fixed amount of melatonin conjugated to horse radish peroxidase (Melatonin-HRP) for binding sites with a rabbit melatonin monoclonal 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 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 Melatonin-HRP detected is inversely proportional to the amount of melatonin in a sample.

## REAGENTS PROVIDED AND REAGENT PREPARATION

Store all other reagents at 2 to 8°C. Use only reagents supplied with this kit. Do not interchange reagents with different lot numbers. Expiration dates and lot numbers are printed on the labels.

- GARGG Plate:** One 96 well microplate (12x8 breakable strip wells) coated with goat anti-rabbit gamma globulin placed in a resealable foil bag with desiccant. One (1) 96 well kit is sufficient for 38 duplicate measurements.
- Concentrated Stock Melatonin (synthetic) solution:** 1 bottle 0.200mL.  
Dilute the 6400 pg/mL stock solution 1:100 (1 part 6400 pg/mL+ 99 parts assay diluent) to obtain the highest working calibrator (64 pg/mL) then, dilute serially 1:2 (starting with the 64 pg/mL calibrator) to obtain the following concentrations of **working calibrators:** 32 pg/mL, 16 pg/mL, 8 pg/mL, 4 pg/mL, 2 pg/mL and 1 pg/mL. "0" calibrator is assay diluent.
- Assay diluent:** 1 bottle 20 mL.
- Stock Melatonin Control Concentrate, 3 ng/mL (3000 pg/mL):** 1 bottle 0.200 mL.

### Working Control #2 preparation (example)

Stock Melatonin Control concentrate 3000 (pg/mL)	Assay Diluent	Dilution	Target	Number of EIA wells per 5 mL volume
0.05 mL (50µL)	4.950 mL	1:100	30 pg/mL	100

### Working control #1 Preparation

Melatonin Control #2	Assay Diluent	Dilution	Target	Number of EIA wells per 5 mL volume
0.5 mL (50µL)	4.5 mL	1:10	3 pg/mL	100



Immediately after use, store the unused portions of the **working calibrators** and the **High** and **Low** Controls at 2-8°C. Discard if not used within **7** days of mixing.

5. **Salivary Melatonin EIA rabbit monoclonal Antibody:** 1 bottle, 3 mL. The solution is blue.
6. **Salivary Melatonin-Horseradish Peroxidase (HRP) concentrate.** 1 amber bottle, 0.7 mL. Melatonin derivative is conjugated to horseradish peroxidase. The solution is yellow and light sensitive.
7. **Melatonin-Horseradish Peroxidase (HRP) conjugate buffer, pH 7.4:** 1 bottle, 6 mL. Use only for the preparation of the **Melatonin-HRP working reagent**.

#### **Melatonin-HRP working reagent preparation:**

Determine the amount of **working Melatonin HRP** needed and dilute 1:10 with **conjugate buffer** pH 7.4. For example, mix 0.5 mL of **Melatonin-HRP concentrate** plus 4.5 mL with **conjugate buffer**. This is sufficient for 100 EIA wells. The **Melatonin-HRP working reagent** is light sensitive. Immediately after use, wrap the vial with the unused portion of the **Melatonin-HRP working reagent** with aluminum foil or alternatively, prepare the **Melatonin-HRP working reagent** in an amber vial. Store at 2-8°C. Discard if not used within 7 days of mixing.

8. **Wash solution (10X concentrated) EIA #1:** 1 bottle, 50 mL of phosphate buffered saline, pH 7.4. Prior to use dilute 1:10 with deionized water.
9. **Color Development Reagent EIA #1:** 1 amber plastic bottle, 15 mL of Tetramethylbenzidine (TMB) plus hydrogen peroxide. Light sensitive.
10. **Stopping Solution EIA #1:** 1 bottle of a 15 mL mixture of diluted sulfuric and hydrochloric acid solution.

#### **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 the foil bag has been opened, care should be taken to reseal tightly.
- Opened kits retain activity for 28 days if stored as described above.
- Expiration dates and lot numbers are printed on the labels.

#### **OTHER MATERIALS REQUIRED, BUT NOT PROVIDED**

- Device to dispense accurately 25 µL and 50 µL.
- Multichannel pipettors.
- Microplate or orbital shaker
- Vortex Mixer
- Microplate washer (not required, plates can be washed manually).
- Microplate reader capable of reading 450 nm with 4 parameter data reduction or comparable software.
- Plate Sealers
- Suitable serum or plasma sample collection device



## SAMPLE COLLECTION AND PREPERATION

- **Collection:** Sample collection and processing procedure must be followed.
- Rinse mouth thoroughly with cold water 5 minutes prior to sample collection
- Do not collect samples when oral diseases, inflammation or lesions exist (blood contamination)
- Saliva can be collected in a suitable sampling device. A minimum of 0.5 mL liquid should be collected.
- After collection, refrigerate sample within 30 minutes and freeze at or below -20°C within 4 hours of collection. On day of assay thaw the saliva samples, vortex and centrifuge at 1500x for 15 minutes. Dispense clear sample into appropriate wells.

## 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 times	<b>TBD</b>

## ASSAY PROCEDURE SUMMARY FLOW CHART

Melatonin Calibrators I.D. (pg/mL)	Calibrator, Control, Sample ( $\mu$ L)	HRP Melatonin Working Reagent ( $\mu$ L)	Anti-Melatonin ( $\mu$ L)	Mix. Incubate for 2 hrs. at Room Temperature, shaking.	Diluted 10X Wash Solution ( $\mu$ L)	Wash 3X	Color Development Reagent ( $\mu$ L)	Mix. Incubate 30 min. at room temperature	Stopping Solution ( $\mu$ L)	Mix. Read at 450nm
0	50	50	25		300		125		125	
1	50	50	25		300		125		125	
2	50	50	25		300		125		125	
4	50	50	25		300		125		125	
8	50	50	25		300		125		125	
16	50	50	25		300		125		125	
32	50	50	25		300		125		125	
64	50	50	25		300		125		125	
Control 1	50	50	25		300		125		125	
Control 2	50	50	25		300		125		125	
Sample	50	50	25		300		125		125	



## ASSAY PROCEDURE

1. It is recommended that the **calibrators, controls** and **samples** should be tested in duplicate and the mean value should be used to report the results.
2. To the GARGG microplate dispense **50 µL** of **working Salivary Melatonin EIA calibrators** (0, 1, 2, 4, 8, 16, 32 and 64 pg/mL), **controls**, and **saliva samples**.
3. Add **50 µL** of **Melatonin-HRP Working Reagent** to all wells.
4. Add **25 µL** of **Anti-Melatonin EIA rabbit monoclonal antibody**.
5. Cover microplate with plastic sealer. Incubate by shaking on a microplate orbital shaker set a 500-900 rpm for **2 hrs.** at room temperature.
6. After incubation, decant the contents of the wells. Wash 3 times with **300 µL** of **diluted wash solution**. After the 3rd wash, invert GARGG microplate on an absorbent paper and tap dry.
7. Dispense **125 µL of Color Development Reagent EIA #1** into each well. Shake briefly (manual). Cover microplate with plastic sealer. Incubate for **30 minutes at room temperature**.
8. Dispense **125 µL of Stopping Solution EIA #1** into each microtiter well of the GARGG plate. Shake briefly (manual). Color changes from blue to yellow.
9. Read at 450 nm on a microplate reader within 10 minutes.

Note: If samples exceed the upper end of the measuring range of 64 pg/mL, dilute with zero calibrator and make appropriate concentration correction.

## TYPICAL RESULTS

Typical Calibration Curve (Actual assay)			
Calibrators (pg/mL)	Mean Absorbance	% B/Bo	Value (pg/mL)
0	2.67	100	0
1	2.41	90.3	1
2	2.21	82.8	2
4	1.95	73.0	4
8	1.53	57.3	8
16	1.01	37.8	16
32	0.60	22.5	32
64	0.36	13.5	64
Control I	2.11	79.0	2.8
Control II	0.67	25.1	29.7
Sample I	2.32	86.9	1.5
Sample II	1.48	55.4	8.2
Sample III	1.08	40.4	15.0



## CALCULATION

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)	1 - 64 pg/mL
----------------------------------	--------------

## QUALITY CONTROL

The expected values for the controls are stated on the certificate of analysis, which are included in the kit. The results can only be accepted if the expected values are met. Follow federal, state and local guidelines for testing quality control materials.

## EXPECTED VALUES

Saliva samples from apparently healthy subjects collected in the PM (bedtime), AM (arise), and noon show the following results below:

Time	Subjects	Median (pg / mL)	Range (pg/mL)
PM	27	6.4	1.4 – 24.2
AM	27	6.3	1.6 – 22.6
NOON	27	2.7	0.2 – 10.2

\*It is recommended that each laboratory establishes its own range of normal values

## PERFORMANCE CHARACTERISTICS

### Specificity of the Antiserum

Compounds	% Cross-Reactivity
N-Acetylserotonin	0.38
5-MethoxyTryptophol	<0.001
5-Methoxy-DL-Thyptophan	<0.001
Serotonin Hydrochloride	<0.001
5-Methoxytryptamine	0.15
6-Hydroxy melatonin	<0.001

## SENSITIVITY

### Analytical Sensitivity

The lower limit of sensitivity was determined by interpolating the mean optical density minus 2 SDs of 20 values at the 0 pg/mL level. The minimal concentration of Melatonin that can be distinguished from 0 is 0.62 pg/mL.

## 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 (pg/mL)	Standard deviation (pg/mL)	%CV
Low	20	5.5	0.275	5.0
Medium	20	9.9	0.507	5.1
High	20	27.6	1.330	4.8



### Inter-assay

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

Sample	N	Mean (pg/mL)	Standard deviation (pg/mL)	%CV
Low	12	5.6	0.620	11.1
Medium	12	9.9	0.971	9.8
High	12	26.6	2.194	8.3

### Inter-lot Variation

The inter-lot precision was determined by duplicate measurements of three (3) serum pools and three (3) individual samples, using three (3) different reagent lots.

Samples ID	Lot # 001 mean (pg/mL)	Lot # 002 mean (pg/mL)	Lot #003 mean (pg/mL)	Inter-lot mean (pg/mL)	Inter-lot Std. Dev. (pg/mL)	Inter-lot CV (%)
Sample 1	23.3	25.7	25.8	24.9	1.415	5.7
Sample 2	11.1	12.9	13.5	12.5	1.249	10.0
Sample 3	4.3	5.1	4.6	4.7	0.404	8.7
Pool 1	2.5	2.7	2.7	2.6	0.115	4.4
Pool 2	18.2	20.9	20.1	19.7	1.387	7.0
Pool 3	34.4	35.8	34.2	34.8	0.872	2.5

### DILUTION STUDY

Sample I.D.	Dilution factor	Expected (pg/mL)	Observed (pg/mL)	Recovery (%)
			<b>27.0</b>	
1	1:2	13.500	13.800	102.2
	1:4	6.750	7.200	106.7
	1:8	3.375	3.300	97.8
	1:16	1.688	1.700	100.7
			<b>17.1</b>	
2	1:2	8.550	9.500	111.1
	1:4	4.275	4.500	105.3
	1:8	2.138	2.200	102.9
	1:16	1.069	1.100	102.9
			<b>23.3</b>	
3	1:2	11.650	10.900	93.6
	1:4	5.825	5.700	97.9
	1:8	2.913	2.900	99.6
	1:16	1.456	1.600	109.9
			<b>39.9</b>	
4	1:2	19.950	21.400	107.3
	1:4	9.975	9.900	99.2
	1:8	4.988	5.600	112.3
	1:16	2.494	2.500	100.3



## RECOVERY

Four saliva samples with different levels of endogenous Melatonin were spiked with known quantities of melatonin

Sample	Endogenous (pg/mL)	Added (pg/mL)	Expected (pg/mL)	Observed (pg/mL)	Recovery (%)
1	11.0	20.0	31.0	33.4	107.7
2	8.0	10.0	18.0	17.0	94.4
3	2.1	20.0	22.1	22.9	103.6
4	17.2	10.0	27.2	29.1	107.0

## LIMITATIONS

- The Melatonin Saliva ELISA Assay Kit reagents are optimized to measure melatonin in human saliva
- Avoid the use of samples containing blood contamination.,
- Samples containing Azide or thimerosal are unsuitable for this assay.
- Avoid repeated freezing and thawing of serum or plasma samples after the initial freeze/thaw.

## PRECAUTIONS

- Only physician, clinical labs, research labs and hospital labs may acquire, possess and use the kit.
- Compare contents and packing list, if there is breakage or shortage, notify Eagle Biosciences immediately.
- Do not pipet reagents by mouth.
- Do not smoke, eat or drink while performing assay.
- Wear disposable rubber gloves.
- Treat all serum or plasma samples as potentially infectious.
- Do not mix reagent lot numbers or alter in any way the reagents in this kit. If this is done, Eagle Biosciences will not be responsible for the performance of the assay.
- Avoid contact with the Color Development Reagent (TMB). It contains solvents that can irritate skin and mucus membranes. If contact is made, wash thoroughly with water.
- Avoid contact with the stopping solution. It contains acid. If contact is made, rinse thoroughly with water.



## REFERENCES

1. Voultsios, A., Kennaway, D.J., & Dawson, D. (1997). Salivary melatonin as a circadian phase marker: Validation and comparison to plasma melatonin. *J. Biol Rhythms*, 12(5), 457-66.
2. Burgess, H.J., & Fogg, L.F. (2008). Individual differences in the amount and timing of salivary melatonin secretion. *PloS One*, 3(8), e3055.
3. Hardeland, R., & Pandi-Perumal, S.R. (2005). Melatonin, a potent agent in antioxidative defense: Actions as a natural food constituent, gastrointestinal factor, drug and prodrug. *Nutr Metab (Lond.)*, 2, 22-36.
4. Tan, D.X., Manchester, L.C., Terron, M.P., Flores, L.J., & Reiter, R. J. (2007). One molecule, many derivatives: A never ending interaction of melatonin with reactive oxygen and nitrogen species? *J. Pineal Res*, 42(1), 28-42.
5. Nowak, R., McMillen, I.C., Redman, J., & Short, R.V. (1987). The correlation between serum and salivary melatonin concentrations and urinary 6hydroxymelatonin sulphate excretion rates: Two non-invasive techniques for monitoring human circadian rhythmicity. *Clin Endocrinol (Oxf.)*, 27(4) 445-
6. deAlmeida, E.A., Di Mascio, P., Harumi, T., Spence, D.W., Moscovitch, A., Hardeland, R., Pandi-Perumal, S.R.(2011). Measurement of melatonin in body fluids: Standards, protocols and procedures. *Childs Nerv Syst*, 27 (6) 879-91

## 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](mailto:info@eaglebio.com) or at 866-411-8023.*