



Norepinephrine ELISA Kit

Catalog Number KA1891

96 assays

Version: 09

Intended for research use only

www.abnova.com

Table of Contents

| | |
|---|-----------|
| Introduction | 3 |
| Intended Use | 3 |
| Principle of the Assay | 3 |
| General Information | 4 |
| Materials Supplied | 4 |
| Storage Instruction | 5 |
| Materials Required but Not Supplied | 5 |
| Precautions for Use | 6 |
| Assay Protocol | 8 |
| Reagent Preparation | 8 |
| Sample Preparation | 8 |
| Assay Procedure | 8 |
| Data Analysis..... | 11 |
| Calculation of Results | 11 |
| Performance Characteristics | 12 |
| Resources | 14 |
| References | 14 |
| Plate Layout | 15 |

Introduction

Intended Use

Enzyme Immunoassay for the quantitative determination of Noradrenaline (Norepinephrine) in plasma and urine.

Principle of the Assay

Noradrenaline (norepinephrine) is extracted by using a cis-diol-specific affinity gel, acylated and then converted enzymatically.

The competitive ELISA kit uses the microtiter plate format. The antigen is bound to the solid phase of the microtiter plate. The derivatized standards, controls and samples and the solid phase bound analytes compete for a fixed number of antibody binding sites. After the system is in equilibrium, free antigen and free antigen-antibody complexes are removed by washing. The antibody bound to the solid phase is detected by an anti-rabbit IgG-peroxidase conjugate using TMB as a substrate. The reaction is monitored at 450 nm. Quantification of unknown samples is achieved by comparing their absorbance with a reference curve prepared with known standard concentrations.

General Information

Materials Supplied

List of component

| Component | Description | Amount |
|---------------------------------|--|-----------------|
| Adhesive Foil | Ready to use. Adhesive foils in a resealable pouch. | 4 slices |
| Wash Buffer Concentrate (50X) | Concentrated buffer with a non-ionic detergent and physiological pH. | 20 mL |
| Enzyme Conjugate | Ready to use, goat anti-rabbit immunoglobulins, conjugated with peroxidase. | 12 mL |
| Substrate | Ready to use, chromogenic substrate containing tetramethylbenzidine, substrate buffer and hydrogen peroxide. | 12 mL |
| Stop Solution | Ready to use, 0.25 M H ₂ SO ₄ . | 12 mL |
| Noradrenaline Microtiter Strips | Ready to use, antigen precoated microwell plate in a resealable yellow pouch with desiccant. | 96 (12x8) wells |
| Noradrenaline Antiserum | Ready to use, rabbit anti-noradrenaline antibody, yellow coloured. | 6 mL |
| Adjustment Buffer | Ready to use. TRIS buffer. | 4 mL |
| Acylation Buffer | Ready to use, buffer with light alkaline pH for the acylation. | 20 mL |
| Acylation Reagent | Ready to use, Acylation reagent in DMF and DMSO. Highly flammable liquid and vapour. May damage fertility or the unborn child. Causes serious eye irritation. | 3 mL |
| Assay Buffer | Ready to use, contains 1 M hydrochloric acid and a non-mercury preservative. | 6 mL |
| Coenzyme | Ready to use, S-adenosyl-L-methionine. | 4 mL |
| Enzyme | Lyophilized, Catechol-O-methyltransferase. | 2 vials |
| Extraction Buffer | Ready to use, buffer containing carbonate. | 6 mL |
| Extraction Plate | Ready to use, coated with boronate affinity gel in a resealable pouch. | 48 wells x 2 |
| Hydrochloric Acid | Ready to use, 0.025 M HCl, yellow coloured. | 20 mL |

Standards and Controls - Ready to use

| Component | Concentration (ng/mL) | Concentration (nmol/mL) | Amount |
|------------|---|-------------------------|--------|
| Standard A | 0 | 0 | 4 mL |
| Standard B | 5 | 30 | 4 mL |
| Standard C | 20 | 118 | 4 mL |
| Standard D | 75 | 443 | 4 mL |
| Standard E | 250 | 1478 | 4 mL |
| Standard F | 1000 | 5910 | 4 mL |
| Control 1 | Refer to QC report for expected value and acceptable range. | | 4 mL |
| Control 2 | Refer to QC report for expected value and acceptable range. | | 4 mL |

Conversion: Noradrenaline (ng/mL) x 5.91 = Noradrenaline (nmol/L)

Content: Acidic buffer with non-mercury stabilizer, spiked with defined quantity of noradrenaline.

Storage Instruction

Store the unopened reagents at 2-8°C until expiration date. Do not use components beyond the expiry date indicated on the kit labels. Once opened the reagents are stable for 1 month when stored at 2-8°C. Once the resealable pouch has been opened, care should be taken to close it tightly with desiccant again.

Materials Required but Not Supplied

- ✓ Calibrated precision pipettes to dispense volumes between 10-700 µL; 1 mL
- ✓ Microtiter plate washing device (manual, semi-automated or automated)
- ✓ ELISA reader capable of reading absorbance at 450 nm and if possible 620-650 nm
- ✓ Microtiter plate shaker (shaking amplitude 3mm; approx. 600 rpm)
- ✓ Absorbent material (paper towel)
- ✓ Water (deionized, distilled, or ultra-pure)
- ✓ Vortex mixer

Precautions for Use

- Procedural cautions, guidelines and warnings
- 1. This kit is intended for professional use only. Users should have a thorough understanding of this protocol for the successful use of this kit. Only the test instruction provided with the kit is valid and has to be used to run the assay. Reliable performance will only be attained by strict and careful adherence to the instructions provided.
- 2. This assay was validated for certain types of samples as indicated in Intended Use. Any off-label use of this kit is in the responsibility of the user and the manufacturer cannot be held liable.
- 3. The principles of Good Laboratory Practice (GLP) have to be followed.
- 4. In order to reduce exposure to potentially harmful substances, wear lab coats, disposable protective gloves and protective glasses where necessary.
- 5. All kit reagents and specimens should be brought to room temperature and mixed gently but thoroughly before use. Avoid repeated freezing and thawing of reagents and specimens.
- 6. For dilution or reconstitution purposes, use deionized, distilled, or ultra-pure water.
- 7. The microplate contains snap-off strips. Unused wells must be stored at 2 °C to 8 °C in the sealed foil pouch with desiccant and used in the frame provided.
- 8. Duplicate determination of sample is highly recommended to be able to identify potential pipetting errors.
- 9. Once the test has been started, all steps should be completed without interruption. Make sure that the required reagents, materials and devices are prepared ready at the appropriate time.
- 10. Incubation times do influence the results. All wells should be handled in the same order and time intervals.
- 11. To avoid cross-contamination of reagents, use new disposable pipette tips for dispensing each reagent, sample, standard and control.
- 12. A standard curve must be established for each run.
- 13. The controls should be included in each run and fall within established confidence limits.
- 14. Do not mix kit components with different lot numbers within a test and do not use reagents beyond expiry date as shown on the kit labels.
- 15. Avoid contact with Stop Solution containing 0.25 M H₂SO₄. It may cause skin irritation and burns. In case of contact with eyes or skin, rinse off immediately with water.
- 16. TMB substrate has an irritant effect on skin and mucosa. In case of possible contact, wash eyes with an abundant volume of water and skin with soap and abundant water. Wash contaminated objects before reusing them.
- 17. For information on hazardous substances included in the kit please refer to Material Safety Data Sheet (MSDS). The Material Safety Data Sheet for this product is made available directly on the website of the manufacturer or upon request.
- 18. The expected reference values reported in this test instruction are only indicative. It is recommended that each laboratory establishes its own reference intervals.

19. Kit reagents must be regarded as hazardous waste and disposed according to national regulations.

- Limitations

Any inappropriate handling of samples or modification of this test might influence the results.

- Interfering substance

- ✓ Plasma: Samples containing precipitates or fibrin strands or which are haemolytic or lipemic might cause inaccurate results.

- ✓ 24-hour urine: please note the sample preparation and storage! If the percentage of the final concentration of acid is too high, the buffer capacity of the Extraction Buffer is insufficient. As a consequence noradrenaline will not be extracted adequately.

- High-Dose-Hook effect

Ni hook effect was observed in this test.

Assay Protocol

Reagent Preparation

- ✓ Wash Buffer: Dilute the 20 mL Wash Buffer Concentrate with water (deionized, distilled, or ultra-pure) to a final volume of 1000 mL.
Storage: 1 month at 2-8°C.

- ✓ Enzyme Solution: Reconstitute the content of the vial labelled 'Enzyme' with 1 mL water (deionized, distilled, or ultra-pure) and mix thoroughly. Add 0.3 mL of Coenzyme followed by 0.7 mL of Adjustment Buffer. The total volume of the Enzyme Solution is 2.0 mL.

Note: The Enzyme Solution has to be prepared freshly prior to the assay (not longer than 10-15 minutes in advance). Discard after use!

Sample Preparation

- Plasma

Whole blood should be collected into centrifuge tubes containing EDTA as anti-coagulant and centrifuged according to manufacturer's instructions immediately after collection. Haemolytic and lipemic samples should not be used for the assay.

Storage: up to 6 hours at 2 - 8°C; for longer periods (up to 6 months) at - 20°C.

Repeated freezing and thawing should be avoided.

- Urine

Spontaneous urine or 24-hour urine, collected in a bottle containing 10-15 mL of 6 M HCl, can be used. If 24-hour urine is used please record the total volume of the collected urine.

Storage: up to 48 hours at 2-8°C, up to 24 hours at room temperature, for longer periods (up to 6 months) at -20°C. Repeated freezing and thawing should be avoided. Avoid exposure to direct sunlight.

Assay Procedure

Allow all reagents to reach room temperature and mix thoroughly by gentle inversion before use. Duplicate determinations are recommended.

The binding of the antiserum and the enzyme conjugate and the activity of the enzyme are temperature dependent, and the absorbance may vary if a thermostat is not used. The higher the temperature, the higher the absorbance will be. Varying incubation times will have a similar influence on the absorbance. The optimal temperature during the Enzyme Immunoassay is between 20-25°C.

- Sample preparation, extraction and acylation
1. Pipette 10 μ L of standards, controls, urine samples and 300 μ L of plasma samples into the respective wells of the Extraction Plate.
 2. Add 250 μ L of water (deionized, distilled, or ultra-pure) to the wells with standards, controls and urine samples.
 3. Pipette 50 μ L of Assay Buffer into all wells.
 4. Pipette 50 μ L of Extraction Buffer into all wells.
 5. Cover plate with adhesive foil and incubate 30 min at RT (20-25°C) on a shaker (approx. 600 rpm).
 6. Remove the foil. Empty plate and blot dry by tapping the inverted plate on absorbent material.
 7. Pipette 1 mL of Wash Buffer into all wells. Incubate the plate for 5 min at RT (20-25°C) on a shaker (approx. 600 rpm). Empty plate and blot dry by tapping the inverted plate on absorbent material.
 8. Pipette another 1 mL of Wash Buffer into all wells. Incubate the plate for 5 min at RT (20-25°C) on a shaker (approx. 600 rpm). Empty plate and blot dry by tapping the inverted plate on absorbent material.
 9. Pipette 150 μ L of Acylation Buffer into all wells.
 10. Pipette 25 μ L of Acylation Reagent into all wells.
 11. Incubate 15 min at RT (20-25°C) on a shaker (approx. 600 rpm).
 12. Empty plate and blot dry by tapping the inverted plate on absorbent material.
 13. Pipette 1 mL of Wash Buffer into all wells. Incubate the plate for 10 min at RT (20-25°C) on a shaker (approx. 600 rpm). Empty plate and blot dry by tapping the inverted plate on absorbent material.
 14. Pipette 150 μ L of Hydrochloric Acid into all wells.
 15. Cover plate with Adhesive Foil. Incubate 10 min at RT (20-25°C) on a shaker (approx. 600 rpm). Remove the foil and discard.

Note: Do not decant the supernatant thereafter!

The following volumes of the supernatant are needed for the subsequent ELISA: Noradrenaline: 20 μ L

- Noradrenaline ELISA
1. Pipette 25 μ L of the Enzyme Solution (refer to Preparation of reagents) into all wells of the Noradrenaline Microtiter Strips.
 2. Pipette 20 μ L of the extracted standards, controls and samples into the appropriate wells.
 3. Incubate for 30 min at RT (20-25°C) on a shaker (approx. 600 rpm).
 4. Pipette 50 μ L of the Noradrenaline Antiserum into all wells and cover plate with Adhesive Foil.
 5. Incubate for 2 hours at RT (20-25°C) on a shaker (approx. 600 rpm).
 6. Remove the foil. Discard or aspirate the content of the wells. Wash the plate 3x by adding 300 μ L of Wash Buffer, discarding the content and blotting dry each time by tapping the inverted plate on absorbent material.
 7. Pipette 100 μ L of the Enzyme Conjugate into all wells.
 8. Incubate for 30 min at RT (20-25°C) on a shaker (approx. 600 rpm).
 9. Discard or aspirate the content of the wells. Wash the plate 3x by adding 300 μ L of Wash Buffer, discarding the content and blotting dry each time by tapping the inverted plate on absorbent material.

10. Pipette 100 μ L of the Substrate into all wells and incubate for 25 ± 5 min at RT (20-25°C) on a shaker (approx. 600 rpm).

Note: Avoid exposure to direct sun light!

11. Add 100 μ L of the Stop Solution to each well and shake the microtiter plate to ensure a homogeneous distribution of the solution.
12. Read the absorbance of the solution in the wells within 10 minutes, using a microplate reader set to 450 nm (if available a reference wavelength between 620 nm and 650 nm is recommended).

Data Analysis

Calculation of Results

- Measuring range of Noradrenaline

Urine: 2.5 to 1000 ng/mL

Plasma: 93-33,333 pg/mL

The standard curve is obtained by plotting the absorbance readings (calculate the mean absorbance) of the standards (linear, y-axis) against the corresponding standard concentrations (logarithmic, x-axis).

Use a non-linear regression for curve fitting (e.g. spline, 4-parameter, akima).

Note: This assay is a competitive assay. This means: the OD-values are decreasing with increasing concentrations of the analyte. OD-values found below the standard curve correspond to high concentrations of the analyte in the sample and have to be reported as being positive.

- Urine samples and controls

The concentrations of the urine samples and the Controls 1 and 2 can be read directly from the standard curve.

Calculate the 24 h excretion for each urine sample: $\mu\text{g}/24\text{h} = \mu\text{g}/\text{L} \times \text{L}/24\text{h}$

- Plasma samples

The read concentrations of the plasma samples have to be divided by 30.

- Conversion

Noradrenaline (ng/mL) \times 5.91 = Noradrenaline (nmol/L)

- Expected reference values

It is strongly recommended that each laboratory should determine its own normal and abnormal values.

24-hour Urine: < 90 $\mu\text{g}/\text{day}$ (535 nmol/day)

Plasma: < 600 pg/mL

- Quality Control

It is recommended to use control samples according to national regulations. Use controls at both normal and abnormal levels. The kit or other commercial controls should fall within established confidence limits.

The confidence limits of the kit controls are printed on the QC-Report.

- Typical standard curve

Example, do not use for calculation!

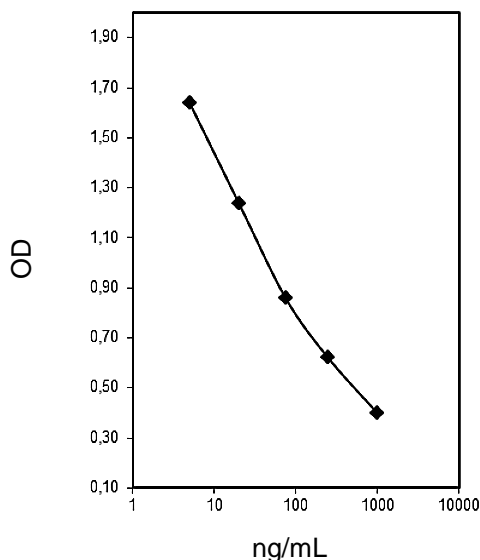


Figure 1: Typical Standard Curve for Norepinephrine ELISA Kit.

Performance Characteristics

- Analytical Sensitivity

| | Sample | Noradrenaline |
|-----|----------------|---------------|
| LOD | Urine (ng/mL) | 1.7 |
| | Plasma (pg/mL) | 0.04 |
| LOQ | Urine (ng/mL) | 2.5 |
| | Plasma (pg/mL) | 93 |

- Analytical Specificity (Cross Reactivity)

| Substance | Cross Reactivity (%) |
|--|----------------------|
| | Noradrenaline |
| Derivatized Adrenaline | 0.08 |
| Derivatized Noradrenaline | 100 |
| Derivatized Dopamine | 0.03 |
| Metanephrine | < 0.01 |
| Normetanephrine | 0.16 |
| 3-Methoxytyramine | < 0.01 |
| 3-Methoxy-4-hydroxyphenylglycol | < 0.01 |
| Tyramine | < 0.01 |
| Phenylalanine, Caffeinic acid, L-Dopa, Homovanillic acid, Tyrosine, 3-Methoxy-4-hydroxymandelic acid | < 0.01 |

- Precision

| Intra-Assay Urine (n=60) | | | |
|--------------------------|--------|---------------|--------|
| | Sample | Range (ng/mL) | CV (%) |
| Noradrenaline | 1 | 26.1 ± 3.6 | 13.8 |
| | 2 | 97 ± 12.8 | 13.4 |
| | 3 | 267 ± 35 | 13.1 |

| Intra-Assay Plasma (n=60) | | | |
|---------------------------|--------|---------------|--------|
| | Sample | Range (pg/mL) | CV (%) |
| Noradrenaline | 1 | 510 ± 65 | 12.8 |
| | 2 | 1358 ± 194 | 14.3 |
| | 3 | 3363 ± 374 | 11.1 |

| Inter-Assay Urine (n=33) | | | |
|--------------------------|--------|---------------|--------|
| | Sample | Range (ng/mL) | CV (%) |
| Noradrenaline | 1 | 19.5 ± 3.9 | 20.0 |
| | 2 | 80.6 ± 10.6 | 13.2 |
| | 3 | 226 ± 39.5 | 17.4 |

| Inter-Assay Plasma (n=18) | | | |
|---------------------------|--------|---------------|--------|
| | Sample | Range (pg/mL) | CV (%) |
| Noradrenaline | 1 | 445 ± 40.9 | 9.2 |
| | 2 | 1232 ± 134 | 10.9 |
| | 3 | 3283 ± 302 | 9.2 |

- Linearity

| | Sample | Serial dilution up to | Range (%) | Mean (%) |
|---------------|--------|-----------------------|-----------|----------|
| Noradrenaline | Urine | 1:512 | 100-127 | 112 |
| | Plasma | 1:512 | 102-125 | 112 |

- Recovery

| | | Mean (%) | Range (%) | Range |
|---------------|--------|----------|-----------|-----------------|
| Noradrenaline | Urine | 103 | 91 – 113 | 58.6-230 ng/mL |
| | Plasma | 87 | 75 – 107 | 51-16,735 pg/mL |

Resources

References

1. Kim et al. Vitamin C prevents stress-induced damage on the heart caused by the death of cardiomyocytes, through the down-regulation of the excessive production of catecholamine, TNF- α , and ROS production in GULO (-/-) Vit C-Insufficient mice. *Free Radical Biology and Medicine*, 65:573-583 (2013).
2. Bada et al. Peripheral vasodilatation determines cardiac output in exercising humans: insight from atrial pacing. *The Journal of Physiology*, 590 (8): 2051-2060 (2012).
3. Parks et al. Employment and work schedule are related to telomere length in women. *Occupational & Environmental Medicine* 68 (8): 582-589 (2011).

Plate Layout

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