

Version 3a, Last updated 15 October 2025

ab211096 Sodium Assay Kit (Colorimetric)

For the sensitive and accurate measurement of sodium in biological fluids.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

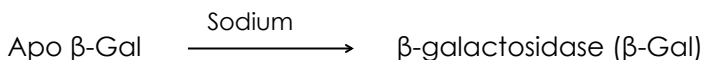
1. Overview	1
2. Protocol Summary	2
3. Precautions	3
4. Storage and Stability	3
5. Limitations	4
6. Materials Supplied	4
7. Materials Required, Not Supplied	4
8. Technical Hints	5
9. Reagent Preparation	6
10. Standard Preparation	7
11. Sample Preparation	8
12. Assay Procedure	9
13. Calculations	10
14. Typical Data	11
15. Quick Assay Procedure	13
16. Troubleshooting	14
17. Interferences	16
18. Notes	17

1. Overview

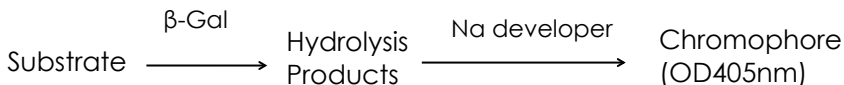
Sodium Assay Kit (Colorimetric) (ab211096) provides a convenient two-step method to detect sodium ions (Na^+) present in serum, urine and saliva. In this assay, sodium ions present in the sample are used by the enzyme β -galactosidase to produce an intermediate product, which reacts with a developer to generate a color signal that can be detected at OD = 405 nm.

This kit can detect as low as 25 μM sodium in biological fluids. Endogenous mono-, di-, and trivalent ions, ascorbic acid, creatinine, glucose, urea, and bilirubin do not interfere with the assay.

Step 1



Step 2



Sodium (Na^+) is one of the most important electrolytes along with chloride, calcium and potassium. Sodium plays vital roles in the maintenance normal cell functions such as plasma volume, pH balance or transmission of nerve impulses. Healthy individuals can absorb sodium ingested in food, and kidneys maintain proper sodium balance by excreting its excess in urine. Normal Sodium intake has been defined to be between 200-500 mg/day.

Hyponatremia (low sodium concentration in blood) can occur in patients with nephrotic syndrome, excessive vomiting and diarrhea, while Hypernatremia (high sodium concentration in blood) is developed in patients suffering from liver diseases, burns, and pregnancy. Traditionally, sodium concentration in clinical settings is determined by potentiometric, gravimetry, photometry, titrimetry and flame atomic emission spectroscopy, but these methods require expensive and complex protocols that need to be performed by trained personnel.

2. Protocol Summary

Standard curve preparation



Sample preparation



Add β -Gal and incubate for 10 minutes at 37°C



Add Substrate and incubate for 30 minutes at 37°C



Measure absorbance (OD405 nm)

3. Precautions

Please read these instructions carefully prior to beginning the assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipette by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

4. Storage and Stability

Store kit at -20°C in the dark immediately upon receipt. Kit has a storage time of 1 year from receipt.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

Aliquot components in working volumes before storing at the recommended temperature.

5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

6. Materials Supplied

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
β -Galactosidase Enzyme	15 μ L	-20°C	-20°C
DTT I (1M)	400 μ L	-20°C	-20°C
Sodium Assay Buffer	25 mL	-20°C	-20°C
Sodium Developer	10 mL	-20°C	-20°C
Sodium Standard	1 mL	-20°C	-20°C
Substrate I	5 mL	-20°C	-20°C

7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring absorbance at OD 405 nm
- MilliQ water or other type of double distilled water (ddH₂O)
- Pipettes and pipette tips, including multi-channel pipette
- Assorted glassware for the preparation of reagents and buffer solutions
- Tubes for the preparation of reagents and buffer solutions
- 96 well plate with clear flat bottom

8. Technical Hints

- **This kit is sold based on number of tests. A “test” simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.**
- Selected components in this kit are supplied in surplus amount to account for additional dilutions, evaporation, or instrumentation settings where higher volumes are required. They should be disposed of in accordance with established safety procedures.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Ensure all reagents and solutions are at the appropriate temperature before starting the assay.
- Samples generating values that are greater than the most concentrated standard should be further diluted in the appropriate sample dilution buffer.
- Make sure all necessary equipment is switched on and set at the appropriate temperature.

9. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.

9.1 β -Galactosidase Enzyme:

Ready to use as supplied. Keep on ice while in use. Aliquot β -Gal so that you have enough volume to perform the desired number of assays. Store at -20°C . Freeze/thaw should be limited. Once opened, use within two months.

9.2 DTT I (1M)

Ready to use as supplied. Equilibrate to room temperature before use. Keep on ice while in use. Store at -20°C .

9.3 Sodium Assay Buffer:

Ready to use as supplied. Equilibrate to room temperature before use. Store at -20°C .

Prior to starting the assay, add DTT I (1M) to Sodium Assay Buffer to a final concentration of 10 mM and label tube as **DTT/Assay Buffer**. Make as much as needed for the assay. Use within 24 hours.

9.4 Sodium Developer:

Ready to use as supplied. Equilibrate to room temperature before use. Store at -20°C .

9.5 Sodium Standard:

Ready to use as supplied. Equilibrate to room temperature before use. Store at -20°C .

9.6 Substrate I:

Ready to use as supplied. Equilibrate to room temperature before use. Mix well before using.

Δ Note: If precipitation is observed in the vial, sonicate the contents in a water bath sonicator at 2-minute interval. Repeat if necessary.

Store at -20°C protected from light. Once opened, use within two months.

10. Standard Preparation

- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.

10.1 Prepare a 7.5 mM Sodium Standard by diluting 5 μL of the 1.5 M Sodium Standard (Step 9.5) in 995 μL of ddH₂O.

10.2 Using 7.5 mM Sodium Standard, prepare standard curve dilutions as described in the table in a microplate or microcentrifuge tubes:

Standard #	Sodium Standard (μL)	DTT/Assay Buffer (μL)	Final volume standard in well (μL)	End amount Sodium in well (nmol/well)
1	0	120	40	0
2	6	114	40	15
3	12	108	40	30
4	18	102	40	45
5	24	96	40	60
6	30	90	40	75

Each dilution has enough amount of standard to set up duplicate readings (2 x 40 μL).

11. Sample Preparation

General sample information:

- We recommend performing several dilutions of your sample to ensure the readings are within the standard value range.
- We recommend that you use fresh samples. If you cannot perform the assay at the same time, we suggest that you snap freeze your samples in liquid nitrogen upon extraction and store them immediately at -80°C. When you are ready to test your samples, thaw them on ice and proceed with the Sample Preparation step. Be aware however that this might affect the stability of your samples and the readings can be lower than expected.
- Sodium concentration range can vary widely. Normal range in humans is 135 – 145 mM for serum, 40 – 220 mmol/day for urine (or > 20 mM for one-time sample), and 30 – 220 mM for saliva.

11.1 Serum:

11.1.1 Dilute serum 1:10 – 1:100 in DTT/Assay Buffer.

Δ Note: Serum samples should not contain any sodium-salt additives such as sodium heparin, sodium EDTA or sodium citrate as they interfere with the results. We recommend using freshly collected serum free of additives or off-the-clot pooled human serum samples.

11.2 Urine:

11.2.1 Dilute urine 1:25 – 1:200 in DTT/Assay Buffer.

11.3 Saliva:

11.3.1 Centrifuge saliva samples at 10,000 x g, 4°C for 10 minutes and collect supernatant.

11.3.2 Dilute supernatant 1:2 – 1:10 in DTT/Assay Buffer.

Δ Note: We suggest using different volumes of sample to ensure readings are within the standard curve range.

12. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
- We recommend that you assay all standards, controls and samples in duplicate.
- Prepare all reagents, working standards, and samples as directed in the previous sections.

Δ Note If you suspect your samples contain substances that can generate background, set up Sample Background Controls.

12.1 Set up reaction wells:

- Standard wells = 40 μ L standard dilutions.
- Sample wells = 1 – 40 μ L samples (adjust volume to 40 μ L/well with DTT/Assay Buffer).
- Background Control Sample wells = 1 – 40 μ L samples (adjust volume to 40 μ L/well with DTT/Assay Buffer).

12.2 β -Gal Reaction:

12.2.1 Dilute β -Galactosidase Enzyme 1:200 by adding 1 μ L of β -Galactosidase Enzyme to 199 μ L of DTT/Assay Buffer. Make as much as needed for the reaction. Keep on ice while in use.

Δ Note: Do not store the diluted β G solution.

12.2.2 Add 20 μ L of diluted β -Galactosidase Enzyme into standard, and sample wells only.

12.2.3 Incubate plate at 37°C for 10 minutes protected from light.

12.2.4 After incubation, add 40 μ L of Substrate I into each well containing Standards, sample background control, and samples. Mix well by pipetting up and down.

12.2.5 Incubate at 37°C for 30 minutes protected from light.

12.2.6 After incubation, add 100 μ L Sodium Developer into each well. Mix well by pipetting up and down.

12.3 Plate measurement:

12.3.1 Measure absorbance on a colorimetric microplate reader at OD = 405 nm in end point mode.

13. Calculations

- Samples producing signals greater than that of the highest standard should be further diluted in appropriate buffer and reanalyzed, then multiply the concentration found by the appropriate dilution factor.

- 13.1** If significant, subtract the sample background control from sample reading.
- 13.2** Average the duplicate reading for each standard and sample.
- 13.3** Subtract the mean absorbance value of the blank (Standard #1) from all standard and sample readings. This is the corrected absorbance.
- 13.4** Plot the corrected absorbance values for each standard as a function of the final concentration of Sodium.
- 13.5** Draw the best smooth curve through these points to construct the standard curve. Most plate reader software or Excel can plot these values and curve fit. Calculate the trendline equation based on your standard curve data (use the equation that provides the most accurate fit).
- 13.6** Apply the corrected sample OD reading to the standard reading to get Sodium (B) amount in the sample wells.
- 13.7** Concentration of Sodium (nmol/ μ L or mM) in the test samples is calculated as:

$$\text{Sodium concentration} = \frac{B}{V} * D$$

Where:

B = amount of Sodium in the sample well calculated from standard curve (nmol).

V = sample volume added in the sample wells (μ L).

D = sample dilution factor.

Sodium Molar Mass = 22.98 g/mol

14. Typical Data

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

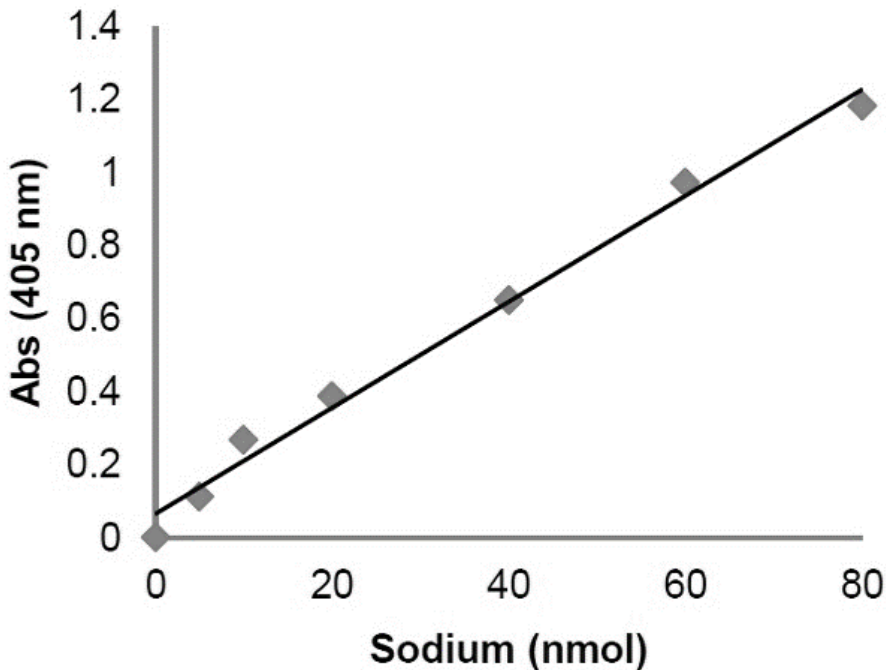


Figure 1. Typical Sodium standard calibration curve.

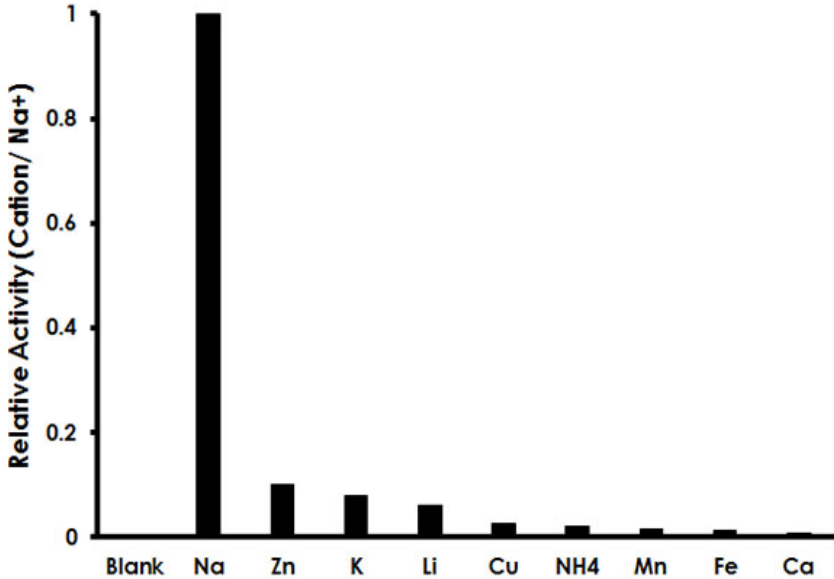


Figure 2. Assay Specificity: Sodium, and other mono, di and trivalent cations (15 nmol/0.15 mM each) were tested to evaluate possible interferences. Interferences were found to be < 10% when data was normalized using Sodium as 100% activity.

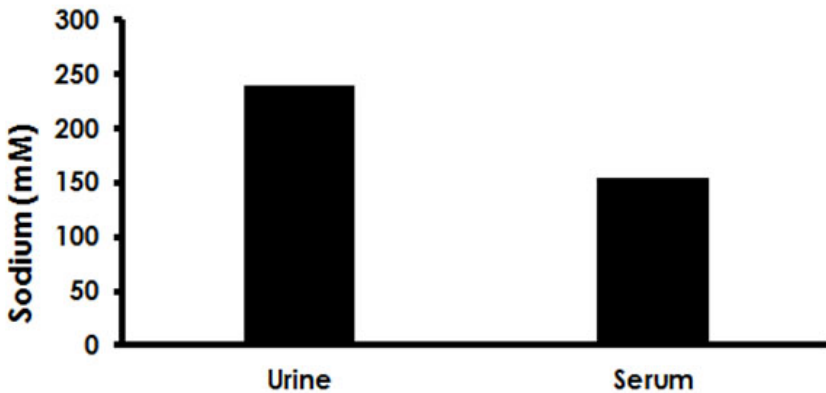


Figure 3. Estimation of sodium in human pooled serum off-the-clot (5 μ L; 1:50 dilution) and human urine (10 μ L; 1:100 dilution).

15. Quick Assay Procedure

Δ Note: this procedure is provided as a quick reference for experienced users. Follow the detailed procedure when performing the assay for the first time.

- Prepare reagents; get equipment ready.
- Prepare DTT/Assay Buffer.
- Prepare Sodium standard dilution [15 – 75 nmol/well].
- Prepare samples in optimal dilutions to fit standard curve readings.
- Set up plate in duplicate for standard (40 μ L), samples (40 μ L) and background sample control wells (40 μ L).
- Dilute β -Galactosidase Enzyme 1:200 in DTT/Assay Buffer.
- Add 20 μ L of diluted β Gal into Standard and sample wells.
- Incubate plate at 37°C for 10 minutes protected from light.
- Add 40 μ L of Substrate I into each well.
- Incubate plate at 37°C for 30 minutes protected from light.
- Add 100 μ L Sodium Developer into each well.
- Measure absorbance on a colorimetric microplate reader at OD = 405 nm in end point mode.

16. Troubleshooting

Problem	Reason	Solution
Assay not working	Use of ice-cold buffer	Buffers must be at assay temperature
	Plate read at incorrect wavelength	Check the wavelength and filter settings of instrument
	Use of a different microplate	Colorimetric: clear plates Fluorometric: black wells/clear bottom plates Luminometric: white wells/clear bottom plates
Sample with erratic readings	Cells/tissue samples not homogenized completely	Use Dounce homogenizer, increase number of strokes
	Samples used after multiple free/ thaw cycles	Aliquot and freeze samples if needed to use multiple times
	Use of old or inappropriately stored samples	Use fresh samples or store at - 80°C (after snap freeze in liquid nitrogen) till use
	Presence of interfering substance in the sample	Check protocol for interfering substances; deproteinize samples
Lower/higher readings in samples and standards	Improperly thawed components	Thaw all components completely and mix gently before use
	Allowing reagents to sit for extended times on ice	Always thaw and prepare fresh reaction mix before use
	Incorrect incubation times or temperatures	Verify correct incubation times and temperatures in protocol

Problem	Reason	Solution
Standard readings do not follow a linear pattern	Pipetting errors in standard or reaction mix	Avoid pipetting small volumes (< 5 μ L) and prepare a master mix whenever possible
	Air bubbles formed in well	Pipette gently against the wall of the tubes
	Standard stock is at incorrect concentration	Always refer to dilutions described in the protocol
Unanticipated results	Measured at incorrect wavelength	Check equipment and filter setting
	Samples contain interfering substances	Troubleshoot if it interferes with the kit
	Sample readings above/ below the linear range	Concentrate/ Dilute sample so it is within the linear range

17. Interferences

These chemical or biological materials will cause interferences in this assay causing compromised results or complete failure:

- Sodium-salt additives such as sodium heparin, sodium EDTA or sodium citrate (for serum collection) as they interfere with the results. We recommend using freshly collected serum free of additives or off-the-clot pooled human serum samples

18. Notes

Technical Support

Copyright © 2025 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)