

Agilent Model 123 Ion Selective Electrode Meter



Introduction

The model 123 ion selective electrode meter is designed for the determination of sodium, potassium, calcium, chloride, and pH in foods, beverages, and other feed and food products. The portable and robust build allows the flexibility of bringing the meter to the samples rather than bringing samples back to the lab. The simple-to-use interface and clear LCD readout allows for quick results with limited training requirements. Make testing quicker, easier, and portable with the Agilent 123 meter.

Protocol: Measuring calcium in juices and other liquids

This procedure is used to measure calcium in fruit juices and other liquids in batches of up to 15 samples for a single calibration and slope.

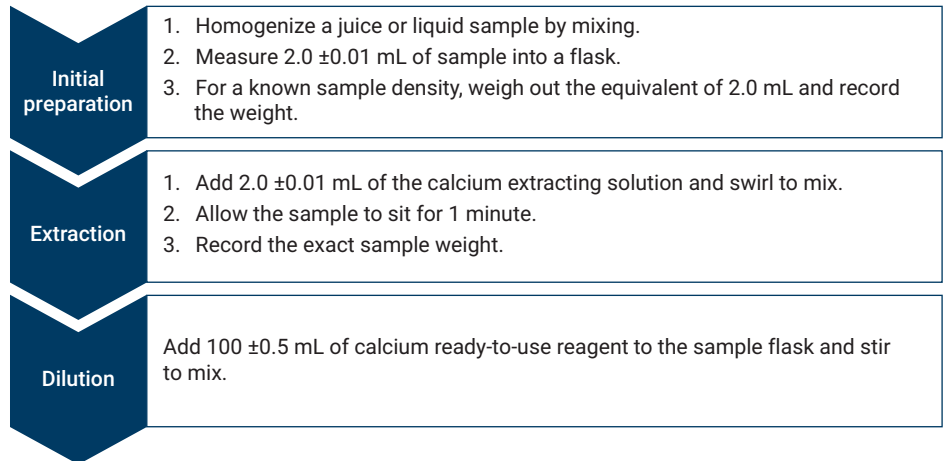
Meter setup

1. Place the meter on a flat surface, away from drafts and out of direct sunlight.
2. Connect the calcium reference electrode (clear plastic barrel) to the electrode jack for Channel 1.
3. Connect the calcium sensing electrode to the electrode BNC connector for Channel 1.
4. Slide the electrodes into the electrode holder and immerse in a sample cup of approximately 40 mL of 100 mg/100 mL calcium calibration standard.
5. Set the *Channel 1 CAL* switch on the back of the meter to 100.
6. Set the *Range* switch to 2000 for Channel 1.
7. Set the *Mode* knob for Channel 1 to Na/K/Ca²⁺.

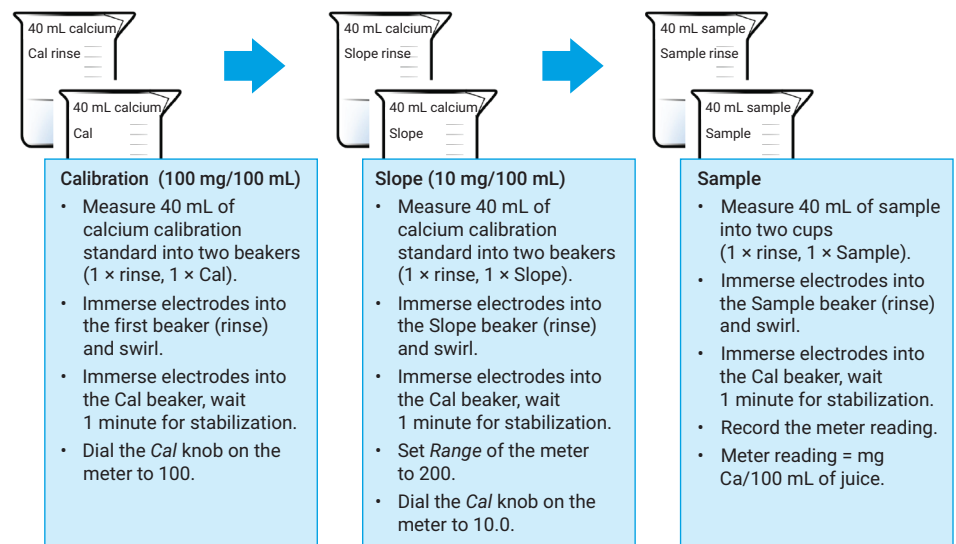
Let the electrodes soak for 12 hours in the standard solution.

| Agilent Part Number | Material Description |
|---------------------|--|
| 123 | Model 123 meter |
| 51770 | Calcium sensing electrode |
| 51010-CA | Calcium reference electrode |
| 82571-G | Calcium ready-to-use reagent |
| 85700-G | Calcium calibration standard, 100 mg/100mL |
| 85790-G | Calcium sloping standard, 10 mg/100mL |
| 55777-Q | Calcium extracting solution |

Sample preparation



Sample analysis



Sample results

| Sample | Volume (mL) | Meter Reading (mg/100 mL) |
|---------------------|-------------|---------------------------|
| Fruit Punch | 2.00 | 4 |
| Apple Juice | 2.00 | 2 |
| Chicken Noodle Soup | 2.00 | 7 |

Protocol: Measuring sodium in liquids and solids

This procedure is used to measure sodium in liquid batches of up to 15 samples for a single calibration and slope.

Meter setup

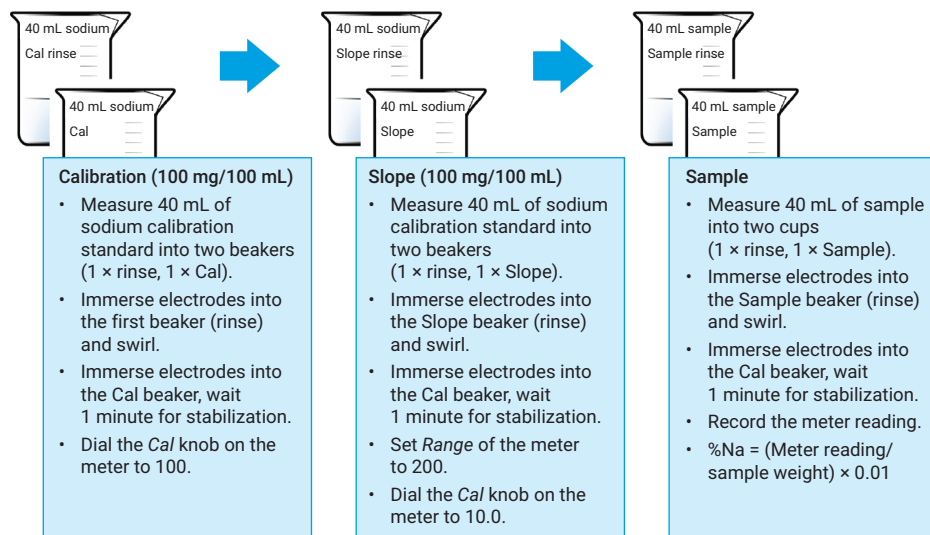
1. Place the meter on a flat surface, away from drafts and out of direct sunlight.
2. Connect the sodium reference electrode (clear plastic barrel) to the electrode jack for Channel 1.
3. Connect the sodium sensing electrode to the electrode BNC connector for Channel 1.
4. Slide the electrodes into the electrode holder and immerse in a sample cup of approximately 40 mL of 100 ppm sodium calibration standard.
5. Set the *Channel 1 CAL* switch on the back of the meter to 100.
6. Set the *Range* switch to 2000 for Channel 1.
7. Set the *Mode* knob for Channel 1 to Na/K/Ca²⁺.

Let the electrodes soak for 12 hours in the standard solution.

Sample preparation

| | |
|---------------------------------------|--|
| Initial liquid preparation | <ol style="list-style-type: none"> 1. Homogenize the juice or liquid sample by mixing. 2. Aliquot 2.0 ±0.01 mL of sample into a flask. 3. For a known sample density, weigh out the equivalent of 2.0 mL and record the weight. |
| Initial solid preparation | <ol style="list-style-type: none"> 1. Homogenize the solid sample with a blender or food processor. 2. Weigh out 1 to 2 g to the nearest 0.01 g of homogenized sample. 3. Record the exact sample weight. |
| Extraction and sample analysis | <ol style="list-style-type: none"> 1. Add 100 ±0.01 mL of the sodium reagent and swirl to mix. 2. Allow the sample to sit for 1 minute. |

Sample analysis



| Agilent Part Number | Material Description |
|---------------------|-------------------------------------|
| 123 | Model 123 Meter |
| 51023 | Sodium sensing electrode |
| 51010-NA | Sodium reference electrode |
| 82506-G | Sodium ready-to-use reagent |
| 82511-G | Sodium calibration standard 100 ppm |
| 82510-G | Sodium sloping standard, 10 ppm |

Sample results

| Sample | Weight/Volume | Meter Reading | % Na |
|---------------------|---------------|---------------|------------|
| Peanut Butter | 5.00 g | 6.5 | 0.01 |
| Fruit Punch | 5.00 mL | 2.5 | Below cal. |
| Apple Juice | 2.00 mL | 6 | 0.03 |
| Chicken Noodle Soup | 2.00 mL | 187 | 0.93 |

Protocol: Measuring chloride in liquids and solids

This procedure is used to measure chloride in liquid batches of up to 15 samples for a single calibration and slope.

Meter setup

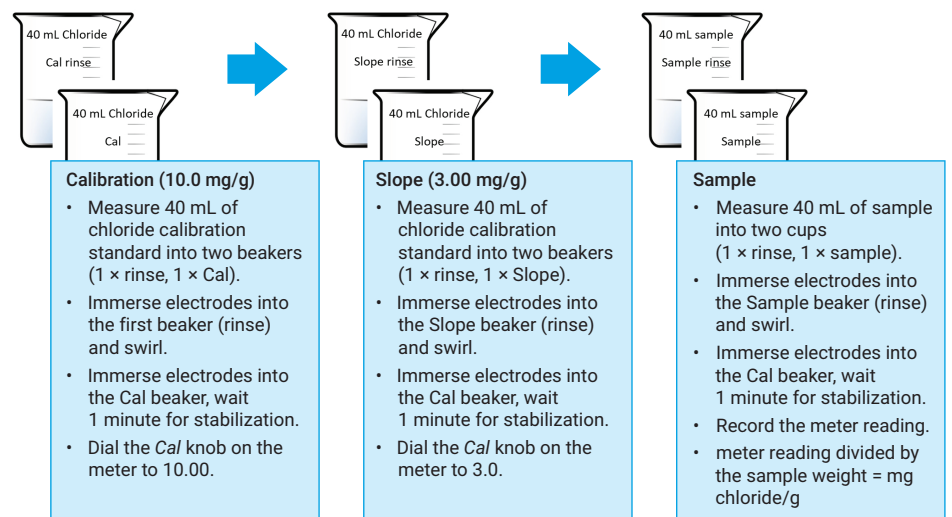
- Place the meter on a flat surface, away from drafts and out of direct sunlight.
- Connect the chloride reference electrode (clear plastic barrel) to the electrode jack for Channel 1.
- Connect the chloride sensing electrode to the electrode BNC connector for Channel 1.
- Slide the electrodes into the electrode holder and immerse in a sample cup of approximately 40 mL of 10 mg/g chloride calibration standard.
- Set the *Channel 1 CAL* switch on the back of the meter to 10.
- Set the *Range* switch to 200 for Channel 1.
- Set the *Mode* knob for Channel 1 to Cl⁻.

Let the electrodes soak overnight in the standard solution.

Sample preparation

| | |
|---------------------------------------|--|
| Initial liquid preparation | <ol style="list-style-type: none"> Homogenize the juice or liquid sample by mixing. Aliquot 2.0 ± 0.01 mL of sample into a flask. For a known sample density, weigh out the equivalent of 2.0 mL and record the weight. |
| Initial solid preparation | <ol style="list-style-type: none"> Homogenize the solid sample with a blender or food processor. Weigh out 1 to 2 g to the nearest 0.01 g of homogenized sample into a beaker. Record the exact sample weight. |
| Extraction and sample analysis | <ol style="list-style-type: none"> Add 100 ± 0.01 mL of the chloride reagent and swirl to mix. Allow the sample to sit for 1 minute. For readings higher than 10, dilute with distilled water. For readings lower than 3, prepare the sample at 5.0 g or 5.0 mL |

Sample analysis



| Agilent Part Number | Material Description |
|---------------------|---------------------------------------|
| 123 | Model 123 meter |
| 51553 | Chloride sensing electrode |
| 51550 | Chloride reference electrode |
| 82551-G | Chloride ready-to-use reagent |
| 82518-G | Chloride calibration standard 10 mg/g |
| 82552-G | Chloride sloping standard, 3 mg/g |

Sample results

| Sample | Volume (mL) | Meter Reading (mg/g) |
|---------------------|----------------------------------|----------------------|
| Fruit Punch | 5.00 | Below calibration |
| Apple Juice | 5.00 | Below calibration |
| Chicken Noodle Soup | 2.00 | 27.3 |
| | 5x dilution with distilled water | 5.5 |

123 Meter parts overview

| Product Number | Description | Category |
|----------------|---|---------------------|
| 123 | Model 123 ISE Meter | Lab-Ready Systems |
| 123-NA | Sodium System | Lab-Ready Systems |
| 123-CL | Chloride System | Lab-Ready Systems |
| 123-K | Potassium System | Lab-Ready Systems |
| 123-CA | Calcium System | Lab-Ready Systems |
| 123-SALT | Salt System | Lab-Ready Systems |
| 123-NA-D | Sodium System with Omnijet Dispenser | Lab-Ready Systems |
| 123-CL-D | Chloride System with Omnijet Dispenser | Lab-Ready Systems |
| 123-K-D | Potassium System with Omnijet Dispenser | Lab-Ready Systems |
| 123-CA-D | Calcium System with Omnijet Dispenser | Lab-Ready Systems |
| 123-SALT-D | Salt System with Omnijet Dispenser | Lab-Ready Systems |
| 54000 | Omnijet Dispenser | Accessories |
| 51023 | Sodium (Na ⁺) Sensing Electrode | Electrodes |
| 51010-NA | Sodium (Na ⁺) Reference Electrode | Electrodes |
| 51553 | Chloride (Cl ⁻) Sensing Electrode | Electrodes |
| 51550 | Chloride (Cl ⁻) Reference Electrode | Electrodes |
| 51333 | Potassium (K ⁺) Sensing Electrode | Electrodes |
| 51010-K | Potassium (K ⁺) Reference Electrode | Electrodes |
| 51770 | Calcium (Ca ⁺) Sensing Electrode | Electrodes |
| 51010-CA | Calcium (Ca ⁺) Reference Electrode | Electrodes |
| 51555 | pH Electrode | Electrodes |
| 82506-G | Sodium/Potassium Reagent RTU – 4 L | Sodium Solutions |
| 82506-5G | Sodium/Potassium Reagent RTU – 20 L | Sodium Solutions |
| 82511-Q | Sodium Standards 100 ppm – 1 L | Sodium Solutions |
| 82511-G | Sodium Standards 100 ppm – 4 L | Sodium Solutions |
| 82511-5G | Sodium Standards 100 ppm – 20 L | Sodium Solutions |
| 82510-Q | Sodium Standards 10 ppm – 1 L | Sodium Solutions |
| 82510-G | Sodium Standards 10 ppm – 4 L | Sodium Solutions |
| 82510-5G | Sodium Standards 10 ppm – 20 L | Sodium Solutions |
| 55555-G | Sodium Extracting Solution – 4 L | Sodium Solutions |
| 55555-5G | Sodium Extracting Solution – 20 L | Sodium Solutions |
| 82507-G | Sodium/Potassium Reagent Concentrate – 4 L | Sodium Solutions |
| 82507-5G | Sodium/Potassium Reagent Concentrate – 20 L | Sodium Solutions |
| 82531-Q | Potassium Standards (100 mg/100 mL) | Potassium Solutions |
| 82532-Q | Potassium Standards (200 mg/100 mL) – 1 L | Potassium Solutions |
| 82532-G | Potassium Standards (200 mg/100 mL) – 4 L | Potassium Solutions |
| 82533-Q | Potassium Standards 50 ppm – 1 L | Potassium Solutions |
| 82533-G | Potassium Standards 50 ppm – 4 L | Potassium Solutions |
| 82555-Q | Potassium Standards 100 ppm – 1 L | Potassium Solutions |
| 82555-G | Potassium Standards 100 ppm – 4 L | Potassium Solutions |

| Product Number | Description | Category |
|----------------|---|--------------------|
| 82557-Q | Salt Standard 1.00 wt% – 1 L | Salt Solutions |
| 82557-G | Salt Standard 1.00 wt% – 4 L | Salt Solutions |
| 82558-Q | Salt Standard 10.00 wt% – 1 L | Salt Solutions |
| 82558-G | Salt Standard 10.00 wt% – 4 L | Salt Solutions |
| 82551-G | Chloride Reagent RTU – 4 L | Chloride Solutions |
| 82551-5G | Chloride Reagent RTU – 20 L | Chloride Solutions |
| 82550-G | Chloride Reagent Concentrate – 4 L | Chloride Solutions |
| 82550-5G | Chloride Reagent Concentrate – 20 L | Chloride Solutions |
| 82518-Q | Chloride Standards (10 mg/g) – 1 L | Chloride Solutions |
| 82518-G | Chloride Standards (10 mg/g) – 4 L | Chloride Solutions |
| 82552-Q | Chloride Standards (3 mg/g) – 1 L | Chloride Solutions |
| 82552-G | Chloride Standards (3 mg/g) – 4 L | Chloride Solutions |
| 82571-G | Calcium Reagent RTU – 4 L | Calcium Solutions |
| 82571-5G | Calcium Reagent RTU – 20 L | Calcium Solutions |
| 82570-G | Calcium Reagent Concentrate – 4 L | Calcium Solutions |
| 82570-5G | Calcium Reagent Concentrate – 20 L | Calcium Solutions |
| 85700-Q | Calcium Standard (100 mg/100 mL) – 1 L | Calcium Solutions |
| 85700-G | Calcium Standard (100 mg/100 mL) – 4 L | Calcium Solutions |
| 85790-Q | Calcium Standard (10 mg/100 mL) – 1 L | Calcium Solutions |
| 85790-G | Calcium Standard (10 mg/100 mL) – 4 L | Calcium Solutions |
| 82533-Q | Calcium Standard 0.10 wt% – 1 L | Calcium Solutions |
| 82573-G | Calcium Standard 0.10 wt% – 4 L | Calcium Solutions |
| 82572-Q | Calcium Standard 1.00 wt% – 1 L | Calcium Solutions |
| 82572-G | Calcium Standard 1.00 wt% – 4 L | Calcium Solutions |
| 82575-Q | Calcium Standard 10.00 wt% – 1 L | Calcium Solutions |
| 82575-G | Calcium Standard 10.00 wt% – 4 L | Calcium Solutions |
| 82572-J-Q | Calcium Standard 1.00 wt% (Fruit Juices) – 1 L | Calcium Solutions |
| 82572-J-G | Calcium Standard 1.00 wt% (Fruit Juices) – 4 L | Calcium Solutions |
| 82575-J-Q | Calcium Standard 10.00 wt% (Fruit Juices) – 1 L | Calcium Solutions |
| 82575-J-G | Calcium Standard 10.00 wt% (Fruit Juices) – 4 L | Calcium Solutions |
| 55777-G | Calcium Extracting Solution – 4 L | Calcium Solutions |
| 55777-5G | Calcium Extracting Solution – 20 L | Calcium Solutions |
| 55400 | Sodium Electrode Etchant Solution | Maintenance |
| 55410 | Sodium/Potassium Reference Fill Solution | Maintenance |
| 55000-P | Potassium Electrode Regenerating Solution | Maintenance |
| 53000-P | Calcium Electrode Regenerating Solution | Maintenance |
| 56000-P | Calcium Reference Fill Solution | Maintenance |
| 52000 | Chloride Electrode Cleaning Compound | Maintenance |
| 55415 | Chloride Reference Fill Solution | Maintenance |
| 55420 | pH Electrode Fill Solution | Maintenance |

Chemical standards portfolio

Save time and order directly online with our new Agilent Chemical Standards Website. Search our catalog of 5,500 environmental, food, and elemental analysis reference materials. Or, order custom standards to meet your specific workflows.

Visit us at www.agilent.com/chem/standards.



www.agilent.com/chem

DE44305.4621412037

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021
Printed in the USA, June 3, 2021
5994-3510EN