

Version 2a Last updated 6 June 2023

ab241008 Nickel Assay Kit

For the measurement of Ni²⁺ in a range of biological samples.

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

PLEASE NOTE: With the acquisition of BioVision by Abcam, we have made some changes to component names and packaging to better align with our global standards as we work towards environmental-friendly and efficient growth. You are receiving the same high-quality products as always, with no changes to specifications or protocols.

Table of Contents

1. Overview	3
2. Materials Supplied and Storage	4
3. Materials Required, Not Supplied	5
4. General guidelines, precautions, and troubleshooting	6
5. Reagent Preparation	7
6. Standard Preparation	8
7. Sample Preparation	9
8. Assay Procedure	10
9. Data Analysis	11
10. Typical Data	13
11. Notes	14

1. Overview

Nickel Assay Kit (ab241008) provides a simple method of quantifying Ni^{2+} in a variety of samples. The assay takes advantage of reaction of Ni^{2+} with mercaptoethanol in borate buffer to form a complex with strong absorbance bands from ~300 to 600 nm. Fe^{2+} and Co^{2+} interfere with the assay, therefore extra steps (as described below) must be taken to subtract the interference in order to determine the correct Nickel concentration in mixed samples. Other ions tested (Mn^{2+} , Cu^{2+} , Zn^{2+}) do not interfere with the assay, presumably no other ionic species are present in high enough concentration to interfere with the reaction. The assay is a simple method of quantifying Ni^{2+} in a variety of samples, which gives a linear range of 2 to 50 nmol nickel containing less than 25 nmol cobalt.

Prepare standards and samples in Nickel Assay Buffer (200 μL final volume).



Add Nickel Reagent (10 μL).



Incubate for 30 minutes at room temperature.



Read OD405 nm (samples without Fe^{2+} or Co^{2+}) or at both OD330 nm and OD405 nm (samples containing Fe^{2+} and/or Co^{2+}).

2. Materials Supplied and Storage

Store kit at room temperature immediately on receipt and check below for storage for individual components. Kit can be stored for 6 months from receipt, if components have not been reconstituted.

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
Nickel Assay Buffer	20 mL	RT	RT
Nickel Reagent	1 mL	RT	RT
Nickel Chloride Standard (1 μ mol)	1 vial	RT	RT

3. 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 330 nm and 505 nm.
- 96 well plate with clear flat bottom

4. General guidelines, precautions, and troubleshooting

Please observe safe laboratory practice and consult the safety datasheet.

For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:

www.abcam.com/assaykitguidelines

For typical data produced using the assay, please see the assay kit datasheet on our website.

5. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.

5.1 Nickel Assay Buffer

Ready to use as supplied.

5.2 Nickel Reagent

Ready to use as supplied.

5.3 Nickel Chloride Standard (1 μmol)

Dissolve in 1 mL dH₂O to make a 1 mM solution. Store at room temperature.

6. Standard Preparation

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

–

- 6.1** Add 0, 10, 20, 30, 40, and 50 μL of 1 mM Nickel Chloride Standard into a series of wells in a 96-well plate to generate 0, 10, 20, 30, 40 and 50 nmol of Nickel Chloride/well. Adjust the volume to 200 μL /well with Nickel Assay Buffer.

Standard #	Nickel Chloride Standard (μL)	Nickel Assay Buffer (μL)	Final volume standard in well (μL)	Nickel Chloride Standard in well (nmol/well)
1	0	200	200	0
2	10	190	200	10
3	20	180	200	20
4	30	170	200	30
5	40	160	200	40
6	50	150	200	50

7. Sample Preparation

General sample information:

We recommend performing several dilutions of your sample to ensure the readings are within the standard value range.

Sample Nickel concentrations can vary over a wide range. Take samples between 10-100 μL and adjust volume to 200 μL with Nickel Assay Buffer for each well. For unknown samples, different sample amounts should be tested to ensure the readings are within the standard curve linear range.

8. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.
- Assay all standards, controls and samples in duplicate.

Δ Note: In the absence of Fe^{2+} and Co^{2+} in samples, the protocol requires reading OD405 nm only. However, in the presence of Fe^{2+} and Co^{2+} in samples, the protocol requires two separate readings at two different wavelengths to correct interference from Fe^{2+} and/or Co^{2+} .

8.1 Reading 1:

Read OD of the samples (200 μL , see Section 7) and standards (200 μL , see Section 6) at 330 nm and 405 nm before adding the Nickel Reagent. This OD is due only to Fe^{2+} and reagent background. Call these measurements OD_{330_1} and OD_{405_1} .

8.2 Development:

Add 10 μL of Nickel Reagent to all standard and sample wells. Incubate at room temperature for 30 minutes to form complex.

8.3 Reading 2:

Read OD at 330 nm and 405 nm. Call these measurements OD_{330_2} and OD_{405_2} .

9. Data Analysis

9.1 Nickel determination in the absence of Fe²⁺ and/or Co²⁺.

Subtract reading 1 (OD405₁) from reading 2 (OD405₂) to get the corrected reading ΔOD405 . Plot the standard curve.

Apply corrected ΔOD405 of unknown samples to the standard curve to determine Ni²⁺ amount in the reaction wells (Ay). Calculate Nickel concentration as in Step 9.3 (without the Step 9.2 corrections).

9.2 Nickel determination in the presence of Fe²⁺ and/or Co²⁺.

Subtract the 0 Nickel OD reading from all standard and sample readings to correct for absorbance due to buffer and plate.

1. **Remove interference at 330 nm due to Fe²⁺** : In the absence of Nickel Reagent, OD330₁ is contributed only by Fe²⁺. After adding Nickel Reagent, Fe²⁺ contribution to the OD330₂ can be calculated as follows: $\text{FeOD330}_2 = 1.82 \times \text{OD330}_1$. Subtract the FeOD330₂ value from total OD330₂, to get corrected OD330, $\Delta\text{FeOD330}_2 = \text{OD330}_2 - \text{FeOD330}_2$, which is contributed by Ni²⁺ and Co²⁺.
2. **Remove interference at 405 nm due to Fe:** In the absence of Nickel Reagent OD405₁ is contributed only by Fe²⁺. After adding Nickel Reagent, Fe²⁺ contribution to OD405₂ can be calculated as follows: $\text{FeOD405}_2 = 1.65 \times \text{OD405}_1$. Subtract FeOD405₂ value from total OD405₂ reading, to get corrected OD405₂ reading, $\Delta\text{FeOD405}_2 = \text{OD405}_2 - \text{FeOD405}_2$, which is contributed by only Ni²⁺ and Co²⁺.
3. **Remove interference due to Co²⁺:** Calculate the ratio of $\Delta\text{FeOD330}_2$ and $\Delta\text{FeOD405}_2$: $\Delta\text{FeOD330}_2/\Delta\text{FeOD405}_2$. The ratio should fall between 0.925 (100% Co) and 2.8125 (100% Ni). Subtract 0.925 from the ratio calculated and divide that result by 1.8875, $((\Delta\text{FeOD330}_2/\Delta\text{FeOD405}_2) - 0.925)/1.8875$ is the percentage of absorbance due to Ni²⁺. Multiply that percentage by $\Delta\text{FeOD330}_2$ to get Nickel absorbance at OD330₂, $\Delta\text{FeCoOD330}_2$ in samples.

9.3 Calculation: Plot the standard curve (ΔOD405 , or ΔOD330).

Calculate sample Nickel reading ΔOD405 from Step 9.1 for samples without Fe²⁺ and Co²⁺, or $\Delta\text{FeCoOD330}_2$ from Step 9.2

for samples with Fe²⁺ or/and Co²⁺ interference. Apply the sample readings to the standard curve to get Ni²⁺ amounts (Ay) in the reaction well.

$$\text{Nickel Concentration (nmol/mL)} = \frac{Ay}{Sv}$$

Where:

Ay = amount of Ni²⁺ in the sample calculated from standard curve (nmol).

Sv = sample volume added in the sample wells (mL).

10. Typical Data

Data provided for demonstration purposes only.

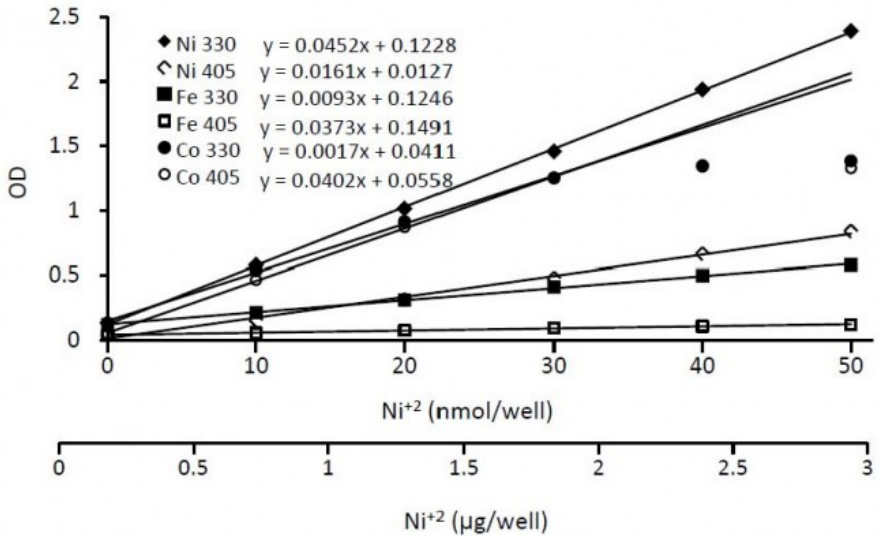


Figure 1. Nickel Standard Curve.

11. Notes

Technical Support

Copyright © 2023 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Austria

wissenschaftlicherdienst@abcam.com | 019-288-259

France

supportscientifique@abcam.com | 01.46.94.62.96

Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

Spain

soportecientifico@abcam.com | 91-114-65-60

Switzerland

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

Canada

ca.technical@abcam.com | 877-749-8807

US and Latin America

us.technical@abcam.com | 888-772-2226

Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

China

cn.technical@abcam.com | +86-21-5110-5938 | 400-628-6880

Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

Singapore

sg.technical@abcam.com | 800 188-5244

Australia

au.technical@abcam.com | +61-(0)3-8652-1450

New Zealand

nz.technical@abc.com | +64-(0)9-909-7829