

# Nickel Colorimetric Assay Kit

(Catalog #K510-100; 100 assays; Store at room temperature)

## I. Introduction:

Nickel is one of four ferromagnetic elements (symbol Ni, at. Num. 28). Several enzymes depend on nickel for activity, including some ureases, carbon monoxide dehydrogenases (methane forming enzymes which reduce CO<sub>2</sub> to CH<sub>4</sub>) and some hydrogenases which allow the production or removal of H<sub>2</sub>. Most of these activities are found in the archaeobacteria. Nickel forms complexes with sulfhydryl compounds with significant absorbance in the UV/visible region in the presence of other ions. BioVision's Nickel Assay kit provides a simple method of quantifying Ni<sup>2+</sup> in a variety of samples. The assay takes advantage of reaction of Ni<sup>2+</sup> with mercaptoethanol in borate buffer to form a complex with strong absorbance bands from ~300 to 600 nm. Fe<sup>2+</sup> and Co<sup>2+</sup> 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<sup>2+</sup>, Cu<sup>2+</sup>, Zn<sup>2+</sup>) 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<sup>2+</sup> in a variety of samples, which gives a linear range of 2 to 50 nmol Nickel containing less than 25 nmol Cobalt.

## II. Kit Contents:

Components	K510-100	Cap Code	Part Number
Nickel Assay Buffer	20 ml	WM	K510 -100-1
Nickel Reagent	1 ml	Green	K510 -100-2
Nickel Chloride Standard (1.0 μmol)	lyophilized	Yellow	K510 -100-3

## III. Reagent Preparation and Storage Conditions:

Store kit at room temperature, keep tightly capped.

**Nickel Assay Buffer and Nickel Reagent:** Ready to use as supplied. Store at room temperature. Stable for at least 6 months.

**Nickel Standard:** Dissolve in 1 ml dH<sub>2</sub>O to make a 1 mM solution. Store at room temperature.

## IV. Nickel Assay Protocol:

- Standard Curve Preparations:** Add 0, 10, 20, 30, 40, 50 μl of the Nickel standard to a series of wells. Adjust volume to 200 μl/well with Nickel Assay Buffer to generate 0, 10, 20, 30, 40 and 50 nmol per well of the Nickel Standard.
- Sample Preparation:** 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.  
**Note:** In the absence of Fe<sup>2+</sup> and Co<sup>2+</sup> in samples, the protocol requires reading OD405 nm only. In the presence of Fe<sup>2+</sup> and Co<sup>2+</sup> in samples, the protocol requires two separate readings at two different wavelengths to correct interference from Fe<sup>2+</sup> and/or Co<sup>2+</sup>.
- Reading 1:** Read OD of the samples and standards at 330 nm and 405 nm before adding the Nickel Reagent. This OD is due only to Fe<sup>2+</sup> and reagent background. Call these measurements OD330<sub>1</sub> and OD405<sub>1</sub>.
- Development:** Add 10 μl of Nickel Reagent to all standard and sample wells. Incubate at room temperature for 30 min to form complex.
- Reading 2:** Read OD at 330 nm and 405 nm. Call these measurements OD330<sub>2</sub> and OD405<sub>2</sub>.
- Nickel Determination in the absence of Fe<sup>2+</sup>, or/and Co<sup>2+</sup>:** Subtract reading 1 (OD405<sub>1</sub>) from reading 2 (OD405<sub>2</sub>) to get the corrected reading ΔOD405. Plot the standard curve. Apply corrected ΔOD405 of unknown samples to the standard curve to determine Ni<sup>2+</sup> amount in the reaction wells (Ay). Calculate Nickel concentration as in step 8 (without the step 7 corrections).

**FOR RESEARCH USE ONLY! Not to be used on humans.**

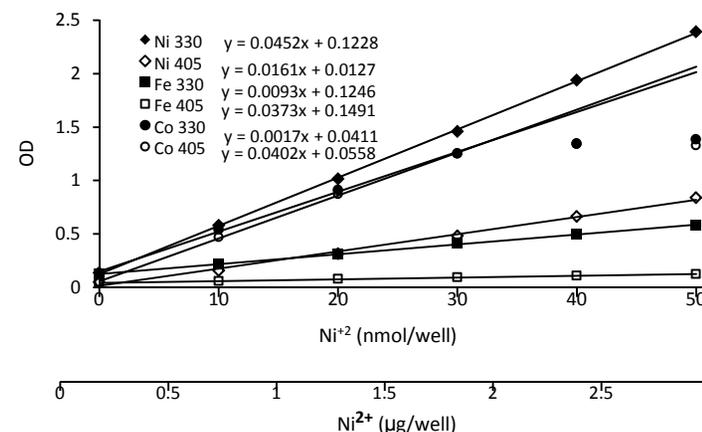
- Nickel Determination in the presence of Fe<sup>2+</sup> or/and Co<sup>2+</sup>:** Subtract the 0 Nickel OD reading from all standard, sample readings to correct absorbance due to buffer and plate.

- Remove interference at 330 nm due to Fe<sup>2+</sup>:** In the absence of Nickel Reagent, OD330<sub>1</sub> is contributed only by Fe<sup>2+</sup>. After adding Nickel Reagent, Fe<sup>2+</sup> contribution to the OD330<sub>2</sub> can be calculated as follows: **FeOD330<sub>2</sub> = 1.82 X OD330<sub>1</sub>**. Subtract the FeOD330<sub>2</sub> value from total OD330<sub>2</sub> to get corrected OD330, ΔFeOD330<sub>2</sub> = OD330<sub>2</sub> – FeOD330<sub>2</sub>, which is contributed by Ni<sup>2+</sup> and Co<sup>2+</sup>.
- Remove interference at 405 nm due to Fe:** In the absence of Nickel Reagent OD405<sub>1</sub> is contributed only by Fe<sup>2+</sup>. After adding Nickel Reagent, Fe<sup>2+</sup> contribution to OD405<sub>2</sub> can be calculated as follows: **FeOD405<sub>2</sub> = 1.65 X OD405<sub>1</sub>**. Subtract FeOD405<sub>2</sub> value from total OD405<sub>2</sub> reading, to get corrected OD405<sub>2</sub> reading, ΔFeOD405<sub>2</sub> = OD405<sub>2</sub> - FeOD405<sub>2</sub>, which is contributed by only Ni<sup>2+</sup> and Co<sup>2+</sup>.
- Remove interference due to Co<sup>2+</sup>:** Calculate the ratio of ΔFeOD330<sub>2</sub> and ΔFeOD405<sub>2</sub>: **ΔFeOD330<sub>2</sub>/ΔFeOD405<sub>2</sub>**. 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, (ΔFeOD330<sub>2</sub>/ΔFeOD405<sub>2</sub> – 0.925)/1.8875 is the percentage of absorbance due to Ni<sup>2+</sup>. Multiply that percentage by ΔFeOD330<sub>2</sub> to get Nickel absorbance at OD330<sub>2</sub>, ΔFeCoOD330<sub>2</sub> in samples.

- Calculation:** Plot the standard curve (ΔOD405, or ΔOD330). Calculate sample Nickel reading ΔOD405 from step 6 for samples without Fe<sup>2+</sup> and Co<sup>2+</sup>, or ΔFeCoOD330<sub>2</sub> from step 7 for samples with Fe<sup>2+</sup> or/and Co<sup>2+</sup> interference. Apply the sample readings to the standard curve to get Ni<sup>2+</sup> amounts (Ay) in the reaction well.

**Nickel Concentration = Ay/Sv (nmol/ml)**

Where: Ay is the amount of Ni<sup>2+</sup> (nmol) in your sample from the standard curve.  
Sv is the sample volume (in ml) added to the sample well.  
Nickel molecular weight: 58.7 g/mol.



**Nickel Standard Curve:** Assays were performed following the kit protocol.

## RELATED PRODUCTS:

- |                         |                       |
|-------------------------|-----------------------|
| Cobalt Assay Kit        | Chloride Assay Kit    |
| Iron Assay Kit          | Phosphate Assay Kit   |
| Calcium Assay Kit       | Magnesium Assay Kit   |
| Ammonia Assay Kit       | Urea Assay Kit        |
| Nitric Oxide Assay Kits | Glutathione Assay Kit |

## GENERAL TROUBLESHOOTING GUIDE:

Problems	Cause	Solution
Assay not working	<ul style="list-style-type: none"> <li>• Use of ice-cold assay buffer</li> <li>• Omission of a step in the protocol</li> <li>• Plate read at incorrect wavelength</li> <li>• Use of a different 96-well plate</li> </ul>	<ul style="list-style-type: none"> <li>• Assay buffer must be at room temperature</li> <li>• Refer and follow the data sheet precisely</li> <li>• Check the wavelength in the data sheet and the filter settings of the instrument</li> <li>• Fluorescence: Black plates (clear bottoms) ; Luminescence: White plates ; Colorimeters: Clear plates</li> </ul>
Samples with erratic readings	<ul style="list-style-type: none"> <li>• Use of an incompatible sample type</li> <li>• Samples prepared in a different buffer</li> <li>• Samples were not deproteinized (if indicated in datasheet)</li> <li>• Cell/ tissue samples were not completely homogenized</li> <li>• Samples used after multiple free-thaw cycles</li> <li>• Presence of interfering substance in the sample (e.g. metal ion chelators)</li> <li>• Use of old or inappropriately stored samples</li> </ul>	<ul style="list-style-type: none"> <li>• Refer data sheet for details about incompatible samples</li> <li>• Use the assay buffer provided in the kit or refer data sheet for instructions</li> <li>• Use the 10 kDa spin cut-off filter or PCA precipitation as indicated</li> <li>• Use Dounce homogenizer (increase the number of strokes); observe for lysis under microscope</li> <li>• Aliquot and freeze samples if needed to use multiple times</li> <li>• Troubleshoot if needed, deproteinize samples</li> <li>• Use fresh samples or store at correct temperatures till use</li> </ul>
Lower/ Higher readings in Samples and Standards	<ul style="list-style-type: none"> <li>• Improperly thawed components</li> <li>• Use of expired kit or improperly stored reagents</li> <li>• Allowing the reagents to sit for extended times on ice</li> <li>• Incorrect incubation times or temperatures</li> <li>• Incorrect volumes used</li> </ul>	<ul style="list-style-type: none"> <li>• Thaw all components completely and mix gently before use</li> <li>• Always check the expiry date and store the components appropriately</li> <li>• Always thaw and prepare fresh reaction mix before use</li> <li>• Refer datasheet &amp; verify correct incubation times and temperatures</li> <li>• Use calibrated pipettes and aliquot correctly</li> </ul>
Readings do not follow a linear pattern for Standard curve	<ul style="list-style-type: none"> <li>• Use of partially thawed components</li> <li>• Pipetting errors in the standard</li> <li>• Pipetting errors in the reaction mix</li> <li>• Air bubbles formed in well</li> <li>• Standard stock is at an incorrect concentration</li> <li>• Calculation errors</li> <li>• Substituting reagents from older kits/ lots</li> </ul>	<ul style="list-style-type: none"> <li>• Thaw and resuspend all components before preparing the reaction mix</li> <li>• Avoid pipetting small volumes</li> <li>• Prepare a master reaction mix whenever possible</li> <li>• Pipette gently against the wall of the tubes</li> <li>• Always refer the dilutions in the data sheet</li> <li>• Recheck calculations after referring the data sheet; Correct for Fe<sup>2+</sup> and Co<sup>2+</sup></li> <li>• Use fresh components from the same kit</li> </ul>
Unanticipated results	<ul style="list-style-type: none"> <li>• Measured at incorrect wavelength</li> <li>• Samples contain interfering substances</li> <li>• Use of incompatible sample type</li> <li>• Sample readings above/below the linear range</li> </ul>	<ul style="list-style-type: none"> <li>• Check the equipment and the filter setting</li> <li>• Troubleshoot if it interferes with the kit</li> <li>• Refer data sheet to check if sample is compatible with the kit or optimization is needed</li> <li>• Concentrate/ Dilute sample so as to be in the linear range</li> </ul>
<p><b>Note:</b> The most probable list of causes is under each problem section. Causes/ Solutions may overlap with other problems.</p>		