

Nickel Assay Kit ab241008

1 Image

Overview

Product name	Nickel Assay Kit
Detection method	Colorimetric
Product 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.
Notes	This product is manufactured by BioVision, an Abcam company and was previously called K510 Nickel Colorimetric Assay Kit. K510-100 is the same size as the 100 test size of ab241008.
Platform	Microplate reader

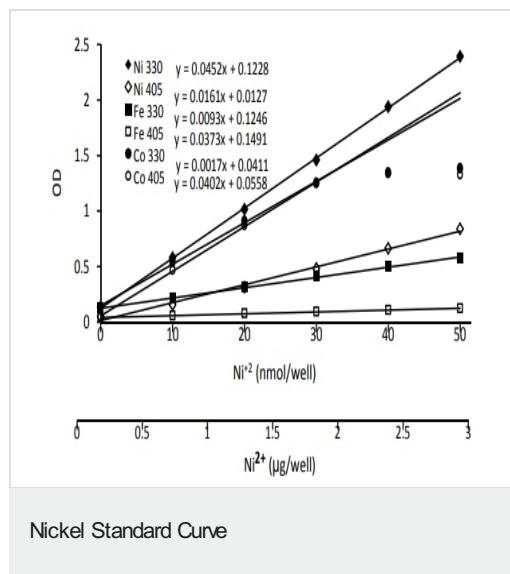
Properties

Storage instructions	Store at room temperature. Please refer to protocols.
----------------------	---

Components	100 tests
Nickel Assay Buffer	1 x 20ml
Nickel Chloride Standard (1.0 µmol)	1 vial
Nickel Reagent	1 x 1ml

Relevance	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 ₂ to CH ₄) and some hydrogenases which allow the production or removal of H ₂ . 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.
-----------	---

Images



Assays were performed following the kit protocol.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors