

Human Nucleoside Diphosphate Kinase D ELISA Kit ab283968

[1 Image](#)

Overview

Product name Human Nucleoside Diphosphate Kinase D ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall	60			5.6%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	60			10.2%

Sample type Cell culture supernatant, Serum, Plasma, Cell Lysate

Assay type Sandwich (quantitative)

Sensitivity 0.34 ng/ml

Range 0.75 ng/ml - 24 ng/ml

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

The Human Nucleoside Diphosphate Kinase D ELISA Kit (ab283968) is designed for detection of Nucleoside Diphosphate Kinase D (NDPKD) in human plasma, serum, cell culture, and cell lysate samples.

This assay employs a quantitative sandwich enzyme immunoassay technique that measures human NDPKD in approximately 5 hours. A polyclonal antibody specific for human NDPKD has been pre-coated onto a 96-well microplate with removable strips. NDPKD in standards and samples is sandwiched by the immobilized antibody and a biotinylated polyclonal antibody specific for human NDPKD, which is recognized by a streptavidin-peroxidase (SP) conjugate. All unbound material is washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions

Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 20ml
1X Standard Diluent	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml
Chromogen Substrate	1 vial
Human Nucleoside Diphosphate Kinase D Microplate	1 unit
Nucleoside Diphosphate Kinase D Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 vial
50X Biotinylated Human Nucleoside Diphosphate Kinase D Antibody	1 vial

Function

Major role in the synthesis of nucleoside triphosphates other than ATP.

Tissue specificity

Widely distributed. Found at very high levels in prostate, heart, liver, small intestine, and skeletal muscle tissues, and in low amounts in the brain and in blood leukocytes.

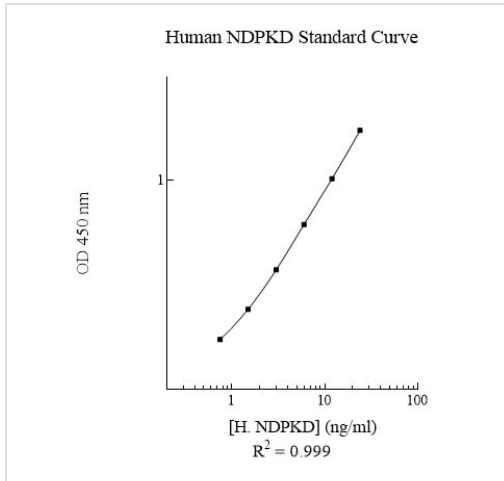
Sequence similarities

Belongs to the NDK family.

Cellular localization

Mitochondrion intermembrane space.

Images



Example of Human Nucleoside Diphosphate Kinase D standard curve. Background-subtracted data values (mean +/- SD) are graphed.

Sandwich ELISA - Human Nucleoside Diphosphate Kinase D ELISA Kit (ab283968)

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