

Human HNRNPK ELISA Kit ab235652

Recombinant SimpleStep ELISA®

5 Images

Overview

Product name Human HNRNPK ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Cell extract	8			7.5%

Sample type

Cell culture extracts

Assay type

Sandwich (quantitative)

Sensitivity

84 pg/ml

Range

117.19 pg/ml - 7500 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture extracts	120	% - %

Assay time

1h 30m

Assay duration

One step assay

Species reactivity

Reacts with: Human

Product overview

Human HNRNPK ELISA Kit (ab235652) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of HNRNPK protein in cell culture extracts. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human HNRNPK with 84 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Notes Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

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

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Human HNRNPK Capture Antibody	1 x 600µl
10X Human HNRNPK Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 4BI	1 x 6ml
Human HNRNPK Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

Function One of the major pre-mRNA-binding proteins. Binds tenaciously to poly(C) sequences. Likely to play a role in the nuclear metabolism of hnRNAs, particularly for pre-mRNAs that contain cytidine-rich sequences. Can also bind poly(C) single-stranded DNA.

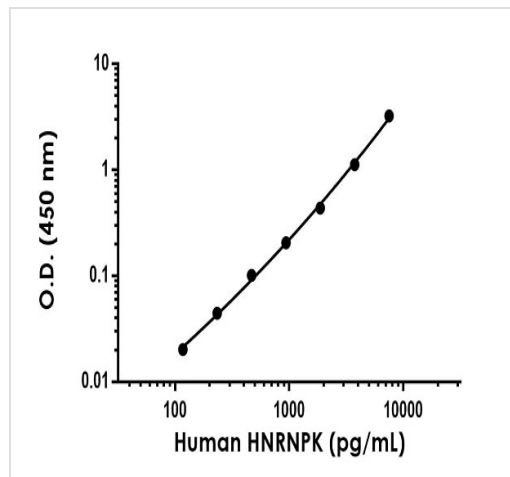
Sequence similarities Contains 3 KH domains.

Post-translational modifications Arg-296 and Arg-299 are dimethylated, probably to asymmetric dimethylarginine.

Cellular localization Cytoplasm. Nucleus > nucleoplasm. In case of ASFV infection, there is a shift in the localization

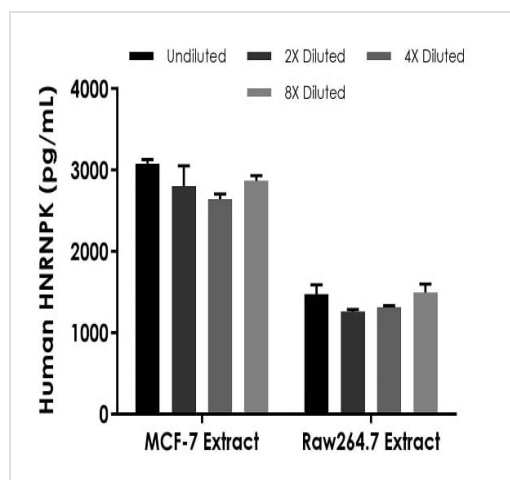
which becomes predominantly nuclear.

Images



Example of human HNRNPK standard curve in 1X Cell Extraction Buffer PTR.

Background-subtracted data values (mean +/- SD) are graphed.



Interpolated concentrations of native HNRNPK in MCF-7 extract and Raw264.7 extract based on a 12.5 $\mu\text{g/mL}$ and 12.5 $\mu\text{g/mL}$ extract loads, respectively.

The concentrations of HNRNPK were measured in duplicate and interpolated from the HNRNPK standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean HNRNPK concentration was determined to be 2,847 pg/mL in MCF-7 extract and 1,384 pg/mL in Raw264.7 extract.

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity

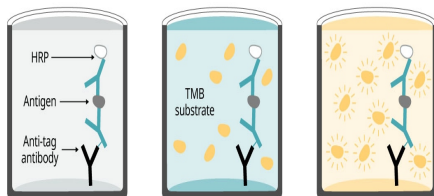


Ethical standards compliant
Animal-free production

Sandwich ELISA - Human HNRNPK ELISA Kit

(ab235652)

To learn more about the advantages of recombinant antibodies see [here](#).



Add sample and antibody cocktail

Wash + add TMB substrate

Add stop solution + Read

90 minutes

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Sandwich ELISA - Human HNRNPK ELISA Kit

(ab235652)

Get more done with SimpleStep ELISA



Easy to use
Single-wash 90-minute protocol



Flexible
Matched antibody pairs available



Precision antibodies
High sensitivity, specificity and reproducibility



Scalable
Now in 10-pack and 384-well formats

Sandwich ELISA - Human HNRNPK ELISA Kit
(ab235652)

To learn more about the advantages of SimpleStep ELISA[®] kits see [here](#).

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