

Product Information & Manual

Hydroxyproline Assay Kit (Colorimetric)
NBP3-24509

Contact

Web: www.bio-technne.com/brands/novus-biologicals/
Email: nb-customerservice@bio-technne.com
P: 888.506.6887 // P: 303.730.1950 // F: 303.730.1966

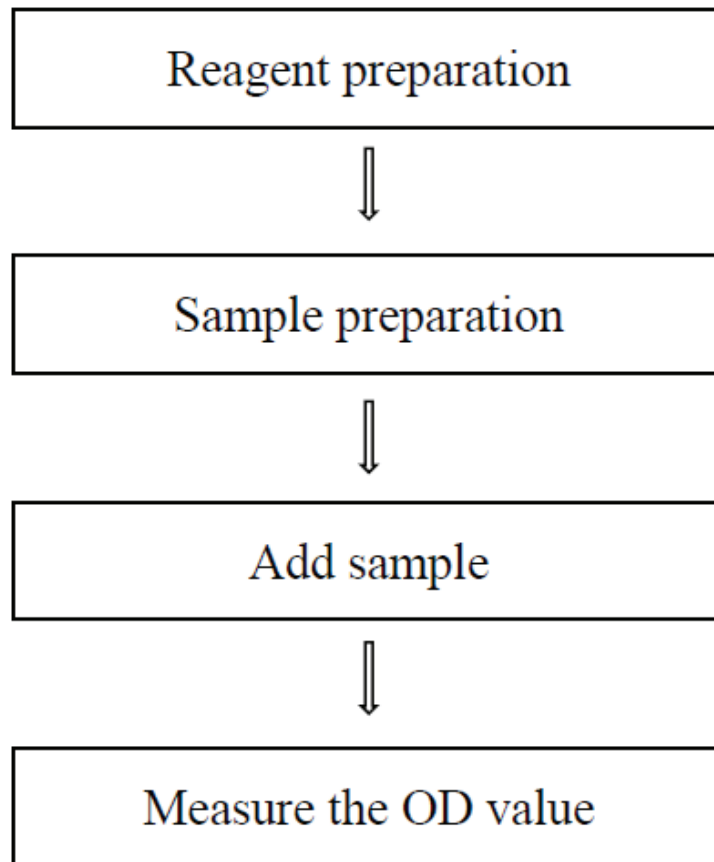
Novus kits are
guaranteed for 6 months
from date of receipt.

**For research use only.
Not for diagnostic or
therapeutic procedures.**

Table of contents

Assay summary	3
Intended use.....	4
Detection principle	4
Kit components & storage	5
Materials prepared by users	5
Reagent preparation	6
Sample preparation.....	7
Operating steps.....	8
Calculation	9
Appendix I Performance Characteristics.....	10
Appendix II Example Analysis	12
Statement.....	13

Assay summary



Intended use

This kit can measure hydroxyproline (HYP) content in serum, plasma and animal tissue samples.

Detection principle

Hydroxyproline is a derivative of proline, mainly found in collagen. Collagen is the most abundant protein in animals and makes up connective tissues such as skin, bone, cartilage, tendons and ligaments. Hydroxyproline enhances the crosslinking of collagen fibers through the hydroxyl group in its structure, thus improving the stability and tensile strength of collagen. Since hydroxyproline almost only exists in collagen, its content can reflect the synthesis and catabolism of collagen. Therefore, the detection of hydroxyproline level can be used as an indicator to monitor collagen metabolism. The detection principle of this kit: the sample is hydrolyzed by the alkali reagent to generate free hydroxyproline, and the oxidation product produced under the action of the oxidant reacts with the color agent to be purple red. The substance has an absorption peak at 565 nm, and within a certain range, its absorbance is linear with the concentration. By measuring the absorbance at 565 nm, the hydroxyproline content in the sample can be calculated.

Kit components & storage

Item	Component	Size 1 (50 Assays)	Size 2 (100 Assays)	Storage
Reagent 1	Buffer Solution	15 mL × 1 vial	30 mL × 1 vial	2-8°C, 12 months, shading light
Reagent 2	Alkali Reagent	50 mL × 1 vial	50 mL × 2 vials	2-8°C, 12 months, shading light
Reagent 3	Oxidant Solution	15 mL × 1 vial	30 mL × 1 vial	2-8°C, 12 months, shading light
Reagent 4	Chromogenic Agent	30 mL × 1 vial	60 mL × 1 vial	2-8°C, 12 months, shading light
Reagent 5	Stop Reagent	30 mL × 1 vial	60 mL × 1 vial	2-8°C, 12 months, shading light
Reagent 6	Clarificant	6 g × 1 vial	6 g × 2 vials	2-8°C, 12 months, shading light
Reagent 7	Acid Reagent	50 mL × 1 vial	50 mL × 2 vials	2-8°C, 12 months, shading light
Reagent 8	100 µg/mL Standard Solution	1.5 mL × 1 vial	1.5 mL × 2 vials	2-8°C, 12 months, shading light

Note: The reagents must be stored strictly according to the preservation conditions in the above table. The reagents in different kits cannot be mixed with each other. For a small volume of reagents, please centrifuge before use, so as not to obtain sufficient amount of reagents.

Materials prepared by users

Instruments:

Spectrophotometer (565 nm), Water bath

Consumptive material:

Precision pH test paper

Reagent preparation

① Equilibrate all reagents to 25°C before use.

② The preparation of measuring working solution:

For each well, prepare 500 μL of measuring working solution (mix well 250 μL of buffer solution and 250 μL of oxidant solution). The measuring working solution should be prepared on spot and used up within the same day.

③ The preparation of 5 $\mu\text{g/mL}$ standard solution:

Before testing, please prepare sufficient 5 $\mu\text{g/mL}$ standard solution. For example, prepare 13 mL of 5 $\mu\text{g/mL}$ standard solution (mix well 0.65 mL of 100 $\mu\text{g/mL}$ standard solution and 12.35 mL of double distilled water). The 5 $\mu\text{g/mL}$ standard solution should be prepared on spot and used up within the same day.

④ The preparation of standard curve:

Always prepare a fresh set of standards. Discard working standard dilutions after use.

Dilute 5 $\mu\text{g/mL}$ standard solution with double distilled water to a serial concentration. The recommended dilution gradient is as follows: 0, 1, 1.5, 2, 3, 3.5, 4, 5 $\mu\text{g/mL}$. Reference is as follows:

Item	①	②	③	④	⑤	⑥	⑦	⑧
Concentration ($\mu\text{g/mL}$)	0	1	1.5	2	3	3.5	4	5
5 $\mu\text{g/mL}$ standard solution (mL)	0	0.6	0.9	1.2	1.8	2.1	2.4	3.0
Double distilled water (mL)	3.0	2.4	2.1	1.8	1.2	0.9	0.6	0

Sample preparation

① Sample preparation

a. Sample hydrolysis:

Tissue sample hydrolysis: accurately weigh 100 mg tissue sample, cut into pieces and put into a glass tube, add 1 mL of alkali reagent, Cover the tube mouth with plastic wrap and open small holes for ventilation, hydrolyze at 95°C for 20 min.

Serum and plasma samples hydrolysis: Add 1 mL of sample and 1 mL of alkali reagent into tube. Cover the tube mouth with plastic wrap and open small holes for ventilation, hydrolyze at 95°C for 20 min.

b. Adjust the pH of the sample hydrolysate:

Cool sample hydrolysate with running water and add 1 mL of acid reagent, mix fully. Add alkali reagent or acid reagent drop by drop. Measure the pH value of the solution to 6.5 using precision pH test paper. Add the double distilled water to a final volume of 10 mL and mix fully.

c. Decolorization of sample hydrolysate:

Take 1.5 mL sample hydrolysate into the centrifugal tube, add about 15 mg of clarificant and mix fully, centrifuge at 15000×g for 10 min, then take the supernatant for detection.

② Dilution of sample

The recommended dilution factor for different samples is as follows (for reference only):

Sample type	Dilution factor
Human serum	1
Rat serum	1
Rat plasma	1
Fish scale	4-10
Fin tissue	4-10
Fishtail tissue	4-10
Mouse lung tissue	1

Mouse muscle tissue	2-5
Mouse leg bone tissue	4-10

Note: The diluent is double distilled water. For the dilution of other sample types, please do pretest to confirm the dilution factor.

Operating steps

- ① Standard tube: Take 1 mL of standard solution with different concentrations to the 5 mL EP tube.
Sample tube: Take 1 mL of sample to the 5 mL EP tube.
- ② Add 0.5 mL of measuring working solution to each tube. Mix fully and stand at room temperature for 20 min.
- ③ Add 0.5 mL of stop reagent to each tube. Mix fully and stand at room temperature for 5 min.
- ④ Add 0.5 mL of chromogenic agent to each tube. Mix fully and water bath at 65°C for 15 min.
- ⑤ Cool the tubes to room temperature with running water. Set the spectrophotometer to zero with double distilled water and measure the OD value of each tube at 565 nm with a 1 cm optical path cuvette.

Calculation

The standard curve:

1. Average the duplicate reading for each standard.
2. Subtract the mean OD value of the blank (Standard #①) from all standard readings. This is the absolved OD value.
3. Plot the standard curve by using absolved OD value of standard and correspondent concentration as y-axis and x-axis respectively. Create the standard curve ($y = ax + b$) with graph software (or EXCEL).

The sample:

1. Serum and plasma samples:

$$\text{HYP content} \begin{matrix} (\mu\text{g/mL}) \end{matrix} = \frac{\Delta A - b}{a} \times V \div V_1 \times f$$

2. Tissue sample:

$$\text{HYP content} \begin{matrix} (\mu\text{g/mg wet weight}) \end{matrix} = \frac{\Delta A - b}{a} \times V \div m \times f$$

[Note]

ΔA : $OD_{\text{Sample}} - OD_{\text{Blank}}$.

V: The volume of sample hydrolysate after pH adjustment, 10 mL.

f: Dilution factor of sample before test.

m: The weight of the sample, mg.

V_1 : The volume of serum (plasma) sample, mL.

Appendix I Performance Characteristics

1. Parameter:

Intra-assay Precision

Three human serum samples were assayed in replicates of 20 to determine precision within an assay (CV = Coefficient of Variation).

Parameters	Sample 1	Sample 2	Sample 3
Mean (µg/mL)	1.00	2.00	4.00
%CV	0.7	1.2	2.0

Inter-assay Precision

Three human serum samples were assayed 20 times in duplicate by three operators to determine precision between assays.

Parameters	Sample 1	Sample 2	Sample 3
Mean (µg/mL)	1.00	2.00	4.00
%CV	3.3	3.5	4.2

Recovery

Take three samples of high concentration, middle concentration and low concentration to test the samples of each concentration for 6 times parallelly to get the average recovery rate of 96%.

	standard 1	standard 2	standard 3
Expected Conc. (µg/mL)	1.00	2.00	4.00
Observed Conc. (µg/mL)	0.95	1.92	3.88
recovery rate(%)	95	96	97

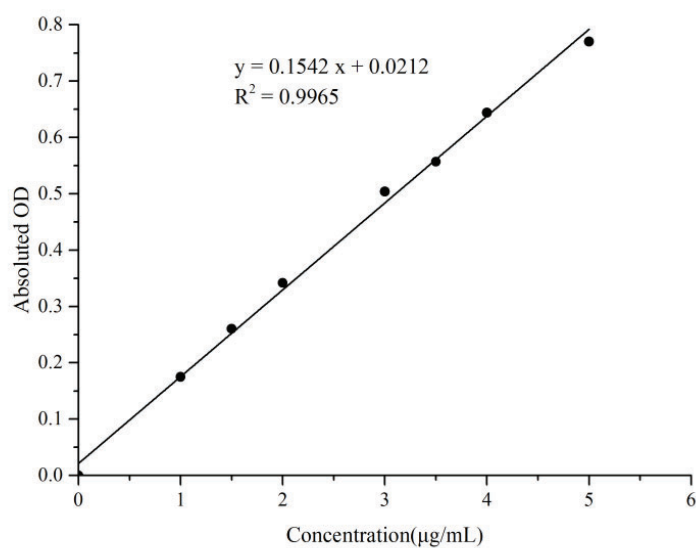
Sensitivity

The analytical sensitivity of the assay is 0.024 µg/mL. This was determined by adding two standard deviations to the mean O.D. obtained when the zero standard was assayed 20 times, and calculating the corresponding concentration.

2. Standard curve:

As the OD value of the standard curve may vary according to the conditions of the actual assay performance (e.g. operator, pipetting technique or temperature effects), so the standard curve and data are provided as below for reference only:

Concentration ($\mu\text{g/mL}$)	0	1	1.5	2	3	3.5	4	5
OD value	0.073	0.248	0.332	0.414	0.578	0.629	0.718	0.844
	0.072	0.247	0.334	0.415	0.575	0.630	0.715	0.841
Average OD value	0.073	0.248	0.333	0.415	0.577	0.630	0.717	0.843
Absoluted OD value	0	0.175	0.261	0.342	0.504	0.557	0.644	0.770



Appendix II Example Analysis

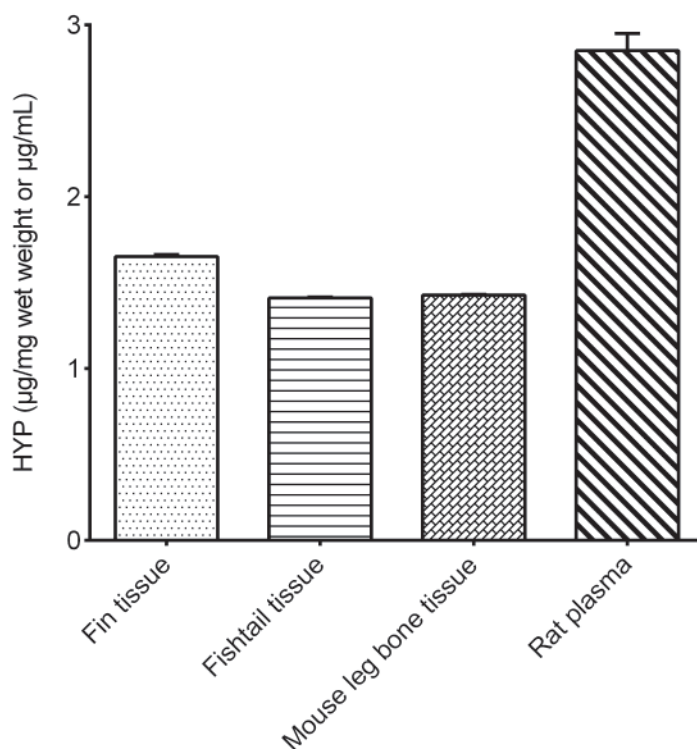
Example analysis:

Take 1 mL of fin tissue which dilute for 4 times and carry the assay according to the operation steps. The results are as follows:

standard curve: $y = 0.1542x + 0.0212$, the average OD value of the blank well is 0.073, the average OD value of the sample well is 0.731, $\Delta A = 0.731 - 0.073 = 0.658$, and the calculation result is:

$$\text{HYP content } (\mu\text{g/mg wet weight}) = \frac{0.658 - 0.0212}{0.1542} \times 10 \times 4 \div 100 = 1.65 \mu\text{g/mg wet weight}$$

Detect fin tissue (The wet weight of sample is 100 mg and the constant volume is 10 mL, dilute for 4 times), fishtail tissue (The wet weight of sample is 100 mg and the constant volume is 10 mL, dilute for 4 times), mouse leg bone tissue (The wet weight of sample is 100 mg and the constant volume is 10 mL, dilute for 4 times) and rat plasma (The volume of sample is 1 mL and the constant volume is 10 mL) according to the protocol, the result is as follows:



Statement

1. This assay kit is for Research Use Only. We will not response for any arising problems or legal responsibilities causing by using the kit for clinical diagnosis or other purpose.
2. Please read the instructions carefully and adjust the instruments before the experiments. Please follow the instructions strictly during the experiments.
3. Protection methods must be taken by wearing lab coat and latex gloves.
4. If the concentration of substance is not within the detection range exactly, an extra dilution or concentration should be taken for the sample.
5. It is recommended to take a pre-test if your sample is not listed in the instruction book.
6. The experimental results are closely related to the situation of reagents, operations, environment and so on. Novus Biologicals will guarantee the quality of the kits only, and NOT be responsible for the sample consumption caused by using the assay kits. It is better to calculate the possible usage of sample and reserve sufficient samples before use.

