

## Product Information

### Hydroxyproline Assay Kit (Perchlorate-Free)

Catalog Number **MAK357**  
Storage Temperature  $-20\text{ }^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

Hydroxyproline ((2*S*,4*R*)-4-hydroxyproline, Hyp) is a nonproteinogenic amino acid formed by the enzyme prolyl-4-hydroxylase, which catalyzes the post-translational modification of polypeptide proline residues following protein synthesis. In animals, hydroxyproline is found almost entirely in collagen, which contains approximately 12–14% hydroxyproline by mass (the fibrillar protein elastin also contains small quantities of hydroxyproline). Hydroxyproline in tissue hydrolysates is thus a direct measure of the amount of collagen present. Collagen is the most abundant protein in mammals and many diseases are believed to affect collagen turnover, including tumor invasion and metastasis, rheumatoid arthritis, cardiopulmonary fibrosis, and muscular dystrophy.

The classical assay for determination of Hyp is based upon oxidation of Hyp to a pyrrole intermediate, followed by reaction with Ehrlich's reagent dissolved in concentrated perchloric acid ( $\text{HClO}_4$ ). Perchloric acid is a hazardous material that is both toxic and highly reactive, requiring special handling and waste-disposal protocols. The Hydroxyproline Assay Kit (Perchlorate-Free) uses a proprietary acidic developer solution to accurately measure Hyp in tissue and protein/peptidase hydrolysates without the use of hazardous perchlorates. It is a quick and convenient protocol, yielding a brightly-colored chromophore with an absorbance peak at 560 nm. The assay can detect a minimum of 0.05  $\mu\text{g}$  of hydroxyproline per well in a 96 well format and generates colorimetric results that are virtually identical to the classical perchloric acid-based method.

This kit is suitable for the estimation of hydroxyproline concentration in tissue homogenates (i.e. muscle, liver, lung, etc.) and human or animal biological fluids (serum, urine).

### Components

The kit is sufficient for 100 colorimetric assays in 96 well plates.

Oxidation Buffer Catalog Number MAK357A	10 mL
Chloramine T Concentrate Catalog Number MAK357B	600 $\mu\text{L}$
Developer Solution Catalog Number MAK357C	5 mL
DMAB Concentrate (in DMSO) Catalog Number MAK357D	5 mL
Hydroxyproline Standard (1 mg/ml) Catalog Number MAK357E	100 $\mu\text{L}$

### Reagents and Equipment Required but Not Provided.

- Sodium Hydroxide (NaOH), 10 M (Catalog Number 221465)
- Hydrochloric Acid (HCl), 37% (12 M) (Catalog Number 258148)
- For hydrolysis: Polypropylene Vials and Screw Caps (Catalog Number TMO362800-0020 or equivalent)
- Pipetting devices and accessories (e.g., multichannel pipettor)
- Centrifuge tubes
- 96 well flat-bottom plate. It is recommended to use clear plates for colorimetric assays
- Spectrophotometric multiwell plate reader

### Precautions and Disclaimer

For R&D use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

### Storage/Stability

The kit is shipped on wet ice. Store components at  $-20\text{ }^{\circ}\text{C}$ , protected from light upon receiving.

- Allow the Oxidation Buffer to warm to room temperature prior to use.
- Chloramine T Concentrate – Provided as a concentrated solution in purified water. Prior to use, warm to room temperature and vortex to ensure it is fully resuspended. Divide into aliquots and store at  $-20\text{ }^{\circ}\text{C}$ , protected from light. Limit aliquots to 3 freeze/thaw cycles.
- Developer Solution – Store at  $4\text{ }^{\circ}\text{C}$  or  $-20\text{ }^{\circ}\text{C}$  and allow solution to come to room temperature prior to use. After use, promptly retighten cap to minimize adsorption of airborne moisture.
- DMAB Concentrate (in DMSO) – Provided as a ready-to-use solution in DMSO. Prior to use, warm to room temperature and mix by vortexing. After use, promptly retighten cap to minimize adsorption of airborne moisture. Store at  $-20\text{ }^{\circ}\text{C}$ , stable for 5 freeze/thaw cycles.
- Hydroxyproline Standard – Provided as a 1 mg/ml stock solution in purified water. Store at  $-20\text{ }^{\circ}\text{C}$ , stable for 5 freeze/thaw cycles.

### Preparation Instructions

#### Reagent Preparation

Oxidation Reagent Mix – For each reaction well to be analyzed (including standard curve wells), prepare 100  $\mu\text{L}$  of Oxidation Reagent Mix by adding 6  $\mu\text{L}$  of Chloramine T Concentrate to 94  $\mu\text{L}$  of Oxidation Buffer. Make a sufficient amount of the Oxidation Reaction Mix for all of the assay wells.

Note: Always prepare fresh Oxidation Reagent Mix as necessary for the number of samples and standards to be quantified. Once diluted to working concentration and exposed to light and air, Chloramine T is only stable for 1–2 hours.

10 M Hydrochloric Acid (HCl) – For every mL of 10 M HCl needed, mix 833  $\mu\text{L}$  of 12 M HCl with 167  $\mu\text{L}$  of ultrapure water.

#### Sample Preparation – Tissue

Add 100  $\mu\text{L}$  of ultrapure water for every 10 mg of tissue and thoroughly homogenize with a glass bead (Dounce) or ultrasonic probe homogenizer. Transfer 100  $\mu\text{L}$  of sample homogenate to a pressure-tight, screw-capped polypropylene vial and add 100  $\mu\text{L}$  of 10 M concentrated NaOH (not provided). Ensure the cap is securely tightened and heat at  $120\text{ }^{\circ}\text{C}$  for 1 hour (See notes). Following alkaline hydrolysis, place vial on ice, allow vial to cool briefly before opening cap and add 100  $\mu\text{L}$  of 10 M concentrated HCl (not provided) to neutralize residual NaOH. Vortex and centrifuge vials at  $10,000 \times g$  for 5 minutes to pellet any insoluble debris that may remain following hydrolysis.

#### Sample Preparation – Urine or Serum

Mix samples with equal volumes of 10 M concentrated NaOH (i.e., 100  $\mu\text{L}$  of urine and 100  $\mu\text{L}$  of 10 M NaOH) in pressure-tight, screw-capped polypropylene vials and hydrolyze at  $120\text{ }^{\circ}\text{C}$  for 1 hour. Cool vials on ice and neutralize hydrolysate by adding an equivalent volume (i.e., 100  $\mu\text{L}$ ) of 10 M concentrated HCl. For urine samples, decolorize with activated charcoal by adding 4 mg of activated charcoal to the neutralized hydrolysate. Vortex and centrifuge at  $10,000 \times g$  for 5 minutes to remove precipitate and activated charcoal, and transfer clarified supernatant to a new tube.

#### Notes:

- Extremely tough samples (containing bone or exoskeletal tissue) may require heating for longer than 1 hour for complete hydrolysis.
- Hydrolysates of certain samples (e.g., fatty tissues) may contain lipid debris that is difficult to pellet by centrifugation. Take care when pipetting hydrolyzed samples to avoid transferring these insoluble globules to sample wells.
- For unknown samples, it is recommended to perform a pilot experiment to ensure readings are within the standard curve range and adjusting the volume of sample hydrolysate accordingly (2–20  $\mu\text{L}$  of hydrolysate may be used) or diluting hydrolysate if necessary.
- For sample hydrolysis, polypropylene vials yield best results.

### Procedure

#### Hydroxyproline Standards

Prepare a 0.1  $\mu\text{g}/\mu\text{L}$  solution of Hydroxyproline Standard by adding 20  $\mu\text{L}$  of the 1 mg/mL Hydroxyproline Standard stock to 180  $\mu\text{L}$  of ultrapure water. Add 0.1  $\mu\text{g}/\mu\text{L}$  solution to a series of wells in a clear, flat-bottom 96 well plate as described in Table 1.

**Table 1.**  
Preparation of Hydroxyproline Standards

Well	0.1 µg/µl Solution	Hydroxyproline (µg/well)
1	0 µL	0
2	2 µL	0.2
3	4 µL	0.4
4	6 µL	0.6
5	8 µL	0.8
6	10 µL	1

#### Assay Reaction

1. Transfer 10 µL of each neutralized sample hydrolysate to desired well(s) in a clear, flat-bottom 96 well plate.
2. Evaporate the sample hydrolysate and standard curve wells to dryness by heating the plate at 65 °C on a hot plate/dry heat block or microplate incubator. To prevent warping/etching of the plastic, do not expose the microplate to extreme temperatures (>85 °C).
3. Add 100 µL of the Oxidation Reagent Mix to each well and incubate the plate at room temperature for 20 minutes.

Note: Following evaporation of the sample hydrolysates, a crystalline residue will be left in the well. Gentle shaking will help dissolve the crystals in Oxidation Reagent Mix more quickly.

Hydrolysates from certain samples may impart a faint yellow tint to the Oxidation Reagent Mix. This slight colorization usually dissipates upon addition of Developer Solution and does not interfere with the assay.

4. Add 50 µL of Developer Solution to each reaction well and incubate the plate at 37 °C for 5 minutes.
  5. Add 50 µL of DMAB Concentrate solution to each reaction well and mix contents thoroughly.
  6. Seal the plate with sealing film and incubate at 65 °C on a hot plate/dry heat block or microplate incubator for 45 minutes.
  7. Measure the absorbance of all sample and standard curve wells at  $A_{560}$  in endpoint mode.
- Note: For maximum signal intensity, measure absorbance within 20 minutes of removing plate from the heat source.

#### Results

For the hydroxyproline standard curve, subtract the zero standard (0 µg/well reagent blank) reading from all standard and test sample readings, plot the background-subtracted values and calculate the slope of the standard curve. For test samples, compare the background-subtracted  $A_{560}$  values to the standard curve to get B µg of hydroxyproline in the well.

Hydrolyzed Hydroxyproline Concentration (µg/µL) =

$$(B/V) \times D$$

where:

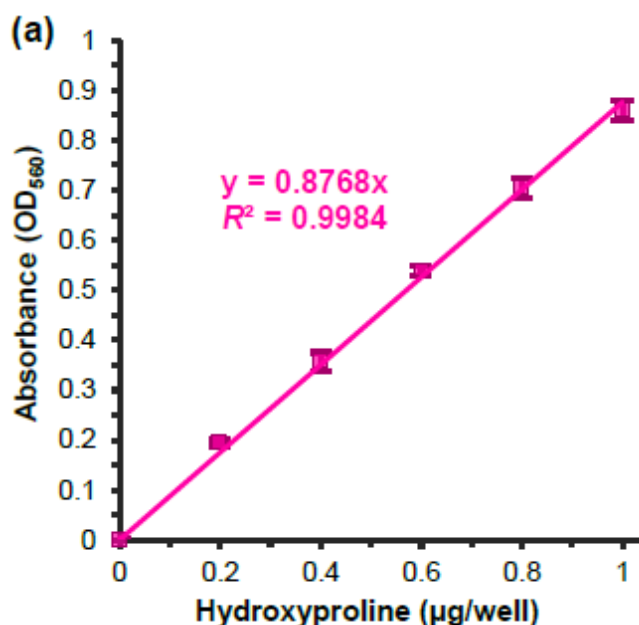
B = The amount of hydroxyproline, calculated from the standard curve (in µg)

V = The volume of sample hydrolysate added to the well (in µL)

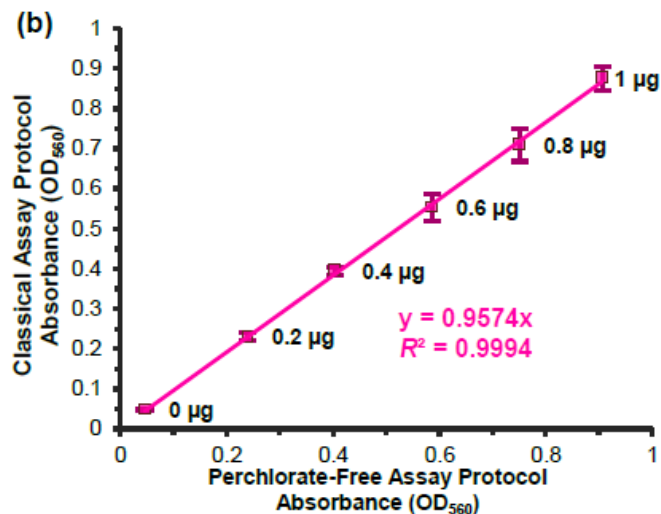
D = The post-hydrolysis sample dilution factor (if applicable, D = 1 for undiluted samples)

Note: The calculation above gives the hydroxyproline concentration in the sample hydrolysate. The dilution factor D is only needed if the sample is diluted after the hydrolysis step. When calculating the amount of hydroxyproline in the original sample homogenate, remember to account for the 3-fold dilution that occurs during generation of the hydrolysate.

**Figure 1.**  
Typical Hydroxyproline Standard Curve

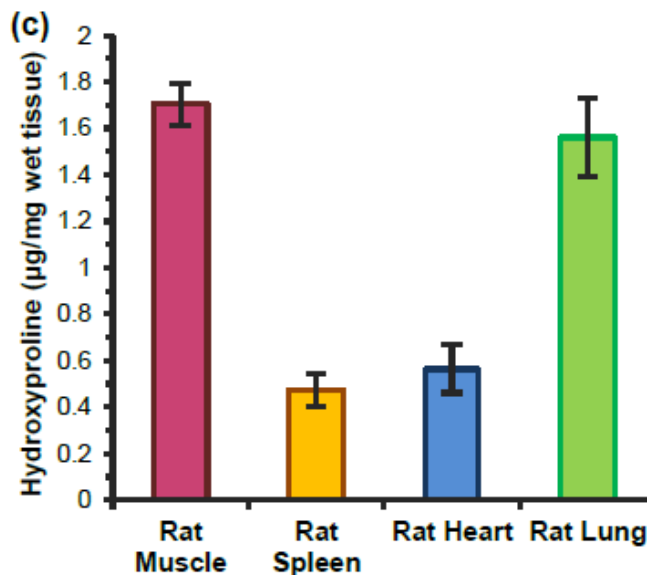


**Figure 2.**  
Comparison of Standard Curves



Correlation between hydroxyproline standard curve absorbance values obtained using the classical perchloric acid-based assay kit and the perchlorate-free assay. The two assay methods show excellent correlation ( $R^2 > 0.999$ ). A Bland-Altman plot indicated a high degree of agreement between the two assays, with a mean difference of only  $3.06 \pm 0.33\%$  between absorbance values generated by the respective methods.

**Figure 3.**  
Hydroxyproline Content in Rat Tissues



Estimation of total hydroxyproline content in rat tissues. Rat leg muscle, spleen, heart, and lung samples were homogenized with ultrapure water, hydrolyzed with 10 M NaOH for 1 hour at 120 °C and neutralized with 10 M HCl. For each sample, 10 µL of the final neutralized hydrolysate was assayed. Hydroxyproline levels (calculated as µg hydroxyproline/mg wet tissue) for the samples were:

muscle,  $1.71 \pm 0.09$  µg/mg  
 spleen,  $0.47 \pm 0.07$  µg/mg  
 heart,  $0.56 \pm 0.10$  µg/mg  
 lung,  $1.56 \pm 0.17$  µg/mg

Data are mean  $\pm$  SEM of 3–4 replicates, assayed according to the kit protocol.

SS, HM, MAM 03/20-1