

Hydroxyproline Assay

A plate-based colorimetric assay to measure 1-20 µg of collagen or about 0.1-2 µg of hydroxyproline in samples.

Hydroxyproline, (MW 131.13), a nonessential amino acid is found primarily in collagen and elastin in animals. Hydroxyproline-rich proteins are also found in the cell walls of plants where it is used as a glycan attachment point. Other instances of hydroxyproline in proteins are known but relatively rare (HIF-1, conotoxin, etc.). Hydroxyproline is formed in an ascorbate dependent process by prolyl hydroxylase after synthesis of the peptide chain. It functions in collagen by strengthening mechanical integrity. Collagen formation and turnover in disease states (liver fibrosis, Paget's, bone metastases) can be assessed through measurement of hydrolysates or serum and urine hydroxyproline levels. AkrivisBio's Hydroxyproline Assay is designed to measure hydroxyproline primarily in protein and tissue hydrolysates. Serum and urine can be used as samples after appropriate pretreatment. AkrivisBio provides a simple method of measuring hydroxyproline giving a chromogen which can be measured between 550-565 nm, useful over the range of 1-20 µg of collagen or about 0.1-2 µg of hydroxyproline.

Ordering Information

[Order Online »](#)

AKR-MA-0101	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

Species Reactivity Species independent

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com