

Product Information Sheet

Product Name: Collagen Type I, FITC-conjugated * Water insoluble*

Catalog Number: AS-85102

Size: 5 mg

Fluorescence: Excitation/Emission wavelength= 490 nm/520 nm

Storage: Collagen type I, FITC-conjugated, is supplied as a lyophilized powder. Store

desiccated at -20°C.

Instruction: Collagen type I is the major collagen existing in skin, bone, tendon and other

fibrous connective tissues. Its fundamental structural unit is tropo-collagen, a 300-kDa triple helix rod formed by three polypeptide α chains. In tendons, tropo-collagens group as collagen fibrils, run parallel to the axis, and have

extensive crosslinkages.

This product is prepared from bovine achilles tendon and then heavily labeled with FITC. It is insoluble in water, aqueous buffers, and organic solvents and may be mechanically minced when preparing it for enzymatic reaction.

A proteolytic degradation assay¹ can be performed by incubation of the FITC labeled collagen type I with collagenase in 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM CaCl₂, for 16-24 h at 37°C. Centrifuge the sample at 10,000 xg for 10 min. Collect the supernatant for fluorescence measurement at Ex/Em=490

nm/520 nm.

For in vitro research use only

References

1. M. N. Sela, D. Kohavi, E. Krausz, D. Steinberg, G. Rosen, *Clin.Oral Implants.Res.* 14, 263-268 (2003).