

ANASPEC

Product Information Sheet

Product Name:	Collagen Type I, FITC-conjugated *Water soluble*
Catalog Number:	AS-85111
Size:	1 mg
Concentration	1mg/mL
Fluorescence:	Excitation/Emission wavelength= 490 nm/520 nm
Storage:	Collagen type I, FITC conjugate, is supplied as 1 mg/mL solution in 0.01 M acetic acid. Store at acid pH at -20°C.
Product description:	<p>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 skin these macromolecules are interlaced and branched to form fibrils. Collagen fibers with limited crosslinkages can partially dissolve in dilute acid.</p> <p>This product is a highly purified telo-peptide free extract from bovine skin, and then heavily labeled with FITC. It is suitable to serve as substrate for collagenases.</p> <p>A proteolytic assay^{1,2} can be performed by incubating the FITC-conjugated collagen type I with collagenases in 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM CaCl₂, at 37°C for 4-16 hr. Centrifuge the reaction solution at 3,000 g for 10 min. Then measure the fluorescence of the supernatant at Ex/Em=490/520 nm.</p> <p>For <i>in vitro</i> research use only</p>

References

1. K. Sugiyama, K. Yamamoto, O. Kamata, N. Katsuda, *Kurume Med J.* 27, 63-69 (1980).
2. A. Baici, G. Cohen, K. Fehr, A. Boni, *Anal.Biochem.* 108, 230-232 (1980).