



Product Information Sheet

Product Name:	Elastin, FITC-conjugated
Catalog Number:	AS-85113
Size:	1 mg
Fluorescence:	Excitation/Emission wavelength= 490 nm/520 nm.
Storage:	Elastin, FITC conjugate, is supplied as lyophilized powder. Store desiccated at -20°C.
Product description:	<p>Elastin soluble is prepared by the method of Partridge.¹ In this method, bovine neck ligament elastin is solublized with hot oxalic acid solution, clarified, dialyzed and lyophilized. The preparation is a salt-free, water soluble powder that readily forms a coacervate at pH 5.0 and 37°C. Elastin is not hydrolyzed by trypsin, chymotrypsin or pepsin but can be cleaved by pancreatic elastase, plant proteases, such as bromelin, ficin and papain and some proteases from microbial sources.</p> <p>Elastin-FITC can be used for elastolytic assays.^{2,3} Incubate the FITC-conjugated elastin with elastases in appropriate assay buffer at 37°C for several hours. The protease activity is demonstrated by the increase of fluorescence in the sample at Ex/Em=490 nm/520 nm.</p> <p>For <i>in vitro</i> research use only</p>

References

1. S.M. Partridge, H.F. Davis, S.T. Adair, *Biochem. J.*, 61, 11 (1955).
2. M. J. Banda, Z. Werb, J. H. McKerrow, *Methods Enzymol.* 144, 288-305 (1987).
3. P. F. Huebner, *Anal.Biochem.* 74, 419-429(1976).