



Product Information Sheet

Product Name: Elastin, FITC-conjugated *Water Insoluble*

Catalog Number: AS-85103

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: Elastin is the insoluble, rubber-like, elastic fibers that are formed by cross-linked polypeptides. It is found in intercellular spaces of the connective tissues of large arteries, trachea, bronchi and ligaments. Elastin is not hydrolyzed by trypsin, chymotrypsin or pepsin but can be cleaved by pancreatic elastase, plant proteases, such as bromelain, ficin and papain and some proteases from microbial sources.

For elastolytic assays^{1,2}, incubate the FITC-conjugated elastin with elastases in appropriate assay buffer at 37°C for 16-24 hr. Centrifuge the sample at 10,000 x g for 10 min. The protease activity is demonstrated by the increase of fluorescence in the supernatant at Ex/Em=490 nm/520 nm.

For *in vitro* research use only

References

1. M. J. Banda, Z. Werb, J. H. McKerrow, *Methods Enzymol.* 144, 288-305 (1987).
2. P. F. Huebner, *Anal. Biochem.* 74, 419-429 (1976).