



Product Information Sheet

Product Name: Gelatin, FITC-conjugated
Catalog Number: AS-85145
Size: 5 mg
Fluorescence: Excitation/Emission wavelength= 495 nm/520 nm.
Storage: Gelatin, FITC conjugate, is supplied as lyophilized powder. Store desiccated at -20°C.

Product description: Gelatin is denatured collagen composed of heterogeneous mixture of water-soluble proteins of high average molecular weights. Below 35-40°C gelatin swells and absorbs 5-10 times its weight of water to form a gel. Gel strength and viscosity gradually weaken upon prolonged heating in solution above 40°C.

This product is extracted from porcine skin, and then heavily labeled with FITC. The stock solution can be prepared by dissolving the product in 10 mM acetic acid with heating to 50°C to facilitate dissolution.

For gelatinolytic assays¹, dilute the stock solution in appropriate assay buffer, pH >7.0. Then incubate the FITC-gelatin with proteases 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 increment of fluorescence in the supernatant at Ex/Em=495 nm/520 nm.

For *in vitro* research use only

References

1. K. Otsuka et al., *J.Nihon Univ Sch Dent.* 39, 182-190 (1997).

ANASPEC INC.

34801 Campus Drive • Fremont, CA 94555 • 510.791.9560 • Fax 510.791.9572
www.anaspec.com