



Product Information Sheet

Product Name:	Collagen Type IV, FAM-conjugated
Catalog Number:	AS-85112
Size:	1 mg
Concentration:	1 mg/mL
Fluorescence:	Excitation/Emission wavelength= 490 nm/520 nm.
Storage:	Collagen type IV, FAM conjugate, is supplied as a solution in 0.01 M acetic acid. Store at acid pH at -20°C.
Product description:	<p>Collagen type IV is the major collagen existing in basement membrane. This product is prepared from human placenta by a modification of the pepsin extraction method¹. The collagen type IV is heavily labeled with 5-FAM and suitable to serve as the substrates for matrix metalloproteinases.</p> <p>For proteolytic assays, dilute the stock solution in appropriate assay buffer, pH >7.0. Then incubate the FAM-collagen type IV 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 increase of fluorescence in the supernatant at Ex/Em=490 nm/520 nm.</p> <p>For <i>in vitro</i> research use only</p>

References

1. Glanville, RW. et al, *Eur.J.Biochem.* 95, 383-389 (1979).