

Casein, FITC-conjugated

Ordering Information

Product Number: 13440 (5 mg)

Storage Conditions

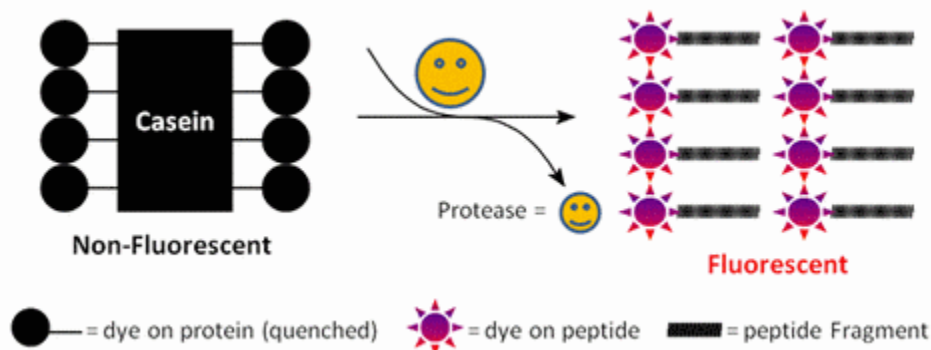
Store at -20 °C

General Properties

Appearance: Light yellow powder
Maximum excitation: 494nm
Maximum Emission: 521 nm
Solvents: Water

Biological Applications

Casein is considered to be a generic substrate for a broad spectrum of proteases. As native casein this fluoresceinated casein is hydrolyzed by many proteases, and widely used for fluorimetric measurement of protease activity. In the intact substrate, casein is heavily labeled with 5-FITC, resulting in significant fluorescence quenching. Protease-catalyzed hydrolysis relieves its quenching effect, yielding brightly green fluorescent dye-labeled short peptides. The increase in fluorescence intensity is directly proportional to protease activity. We do not recommend that this conjugate be used for fluorescence polarization assay. For fluorescence polarization we can custom-make the lightly labeled fluorescein casein conjugate.



Sample Protocol For Trypsin

1. Make a **5-10mg/mL Casein, FITC-conjugated stock solution** in PBS buffer. Unused stock solution can be divided into single use aliquots and stored at -20 °C, and avoid exposure to light.
2. Prepare **2X assay working solution** by diluting the FITC-conjugated stock solution into 50-100 mM Tris buffer (pH 7.4) at 100-400 µg/mL.

Note1: The 2X Assay working solution is designed for detecting the activity of chymotrypsin, trypsin, thermolysin, proteinase K, protease XIV, and human leukocyte elastase. For other proteases, please refer to Appendix I for the appropriate assay buffer formula.

Note2: The optimum concentration of the assay working solution should be determined experimentally for individual proteases.

3. Mix equal volume of the trypsin standards or samples with 2X Assay working solution.
4. Monitor the fluorescence increase at Ex/Em = 490/525 nm.
For kinetic reading: Immediately start measuring fluorescence intensity continuously and record data every 5 minutes for 30 minutes.
For end-point reading: Incubate the reaction at a desired temperature for 30 to 60 minutes, protected from light. Then measure the fluorescence intensity.

Appendix I

<i>Protease</i>	<i>1X Assay Buffer*</i>
Cathepsin D	20 mM Sodium Citrate, pH 3.0
Papain	20 mM sodium acetate, 20 mM cysteine, 2 mM EDTA, pH 6.5
PAE	20 mM sodium phosphate, pH 8.0
Pepsin	10 mM HCl, pH 2.0
Porcine pancreas elastase	10 mM Tris-HCl, pH 8.8
Subtilisin	20 mM potassium phosphate buffer, pH 7.6, 150 mM NaCl

References

1. Vineyard D, Zhang X, Lee I. (2006) Transient kinetic experiments demonstrate the existence of a unique catalytic enzyme form in the peptide-stimulated ATPase mechanism of Escherichia coli Lon protease. *Biochemistry*, 45, 11432.
2. Yadav SC, Pande M, Jagannadham MV. (2006) Highly stable glycosylated serine protease from the medicinal plant Euphorbia milii. *Phytochemistry*, 67, 1414.
3. Frohbieter KA, Ismail B, Nielsen SS, Hayes KD. (2005) Effects of Pseudomonas fluorescens M3/6 bacterial protease on plasmin system and plasminogen activation. *J Dairy Sci*, 88, 3392.
4. Chauhan V, Sheikh AM, Chauhan A, Spivack WD, Fenko MD, Malik MN. (2005) Fibrillar amyloid beta-protein inhibits the activity of high molecular weight brain protease and trypsin. *J Alzheimers Dis*, 7, 37.
5. Lee EH, Kim CS, Cho JB, Ahn KJ, Kim KH. (2003) Measurement of protease activity of live Uronema marinum (Ciliata: Scuticociliatida) by fluorescence polarization. *Dis Aquat Organ*, 54, 85.
6. Cilenti L, Lee Y, Hess S, Srinivasula S, Park KM, Junqueira D, Davis H, Bonventre JV, Alnemri ES, Zervos AS. (2003) Characterization of a novel and specific inhibitor for the pro-apoptotic protease Omi/HtrA2. *J Biol Chem*, 278, 11489.

Disclaimer: This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.