

EnzChek® Peptidase/Protease Assay Kit (E33758)

Quick Facts

Storage upon receipt:

- $\leq -20^{\circ}\text{C}$
- Protect from light
- Avoid freeze/thaw cycles

Ex/Em: 502/528 nm

Introduction

The EnzChek® Peptidase/Protease Assay Kit (E33758) provides a FRET (fluorescence resonance energy transfer)-based method for the simple and accurate quantitation of a wide range of protease activities. The EnzChek® peptidase/protease substrate comprises a fluorophore and a quencher moiety separated by an amino acid sequence. Upon sequence cleavage by protease(s), the fluorophore separates from the quencher and is free to emit a detectable fluorescent signal (excitation and emission maxima of 502 and 528 nm, respectively). The magnitude of the resultant signal is proportional to the degree of substrate cleavage, and can therefore be used to quantitate the enzyme activity present. The assay is performed in a simple mix–incubate–read format, can be completed in 60 minutes or less, and is easily adaptable to accommodate diverse pH requirements.

Materials

Kit Contents

- **Component A:** EnzChek® peptidase/protease substrate, 1.57 mg
- **Component B:** 20X digestion buffer, 3.5 mL of 200 mM Tris-HCl (pH 7.8)
- **Component C:** Substrate solvent, 1 mL of 50% DMSO in 10 mM Tris-HCl (pH 7.8)

Sufficient materials are supplied for 100 assays, based on a 100 μL assay volume in a 96-well microplate format. The EnzChek® peptidase/protease assay can be adapted for use in cuvettes or 384-well microplates.

Storage

Upon receipt, store the kit at $\leq -20^{\circ}\text{C}$, protected from light. Under these conditions the components should be stable for at least 2 years. For short-term storage (days), the buffer (Component B) may be left at room temperature; however, for longer periods we recommend storage at $2-6^{\circ}\text{C}$ to prevent microbial contamination.

Handling and Disposal

We must caution that no data are available addressing the toxicity of the EnzChek® peptidase/protease substrate. Treat the reagent with the same safety precautions as all other chemicals with unknown toxicity, and dispose of the dye in accordance with local regulations.

Protocol

During all steps, protect the EnzChek® peptidase/protease substrate (both concentrate and working solution) from light as much as possible. Allow the kit components to equilibrate to room temperature prior to use.

- 1. Prepare a 1X working solution of digestion buffer .** For example, to prepare enough 1X digestion buffer for ~20 assays, dilute 0.4 mL Component B into 7.6 mL deionized H_2O . This will make 2 mL of 1X digestion buffer available for substrate preparation and enzyme titration, with 6 mL available for intermediate dilutions. Excess buffer is provided for convenient sample preparation. Store the 1X digestion buffer at $2-6^{\circ}\text{C}$.
- 2. Make a stock solution of substrate.** Add 1 mL of substrate solvent (Component C) to the vial containing the EnzChek® peptidase/protease substrate (Component A). Brief sonication may be necessary. For the best results, substrate stock solution should be used within one week of preparation. Unused stock solution may be stored at $\leq -20^{\circ}\text{C}$, but as a result, background signal in the assay may increase over time.
- 3. Prepare a 1X working solution of the EnzChek® peptidase/protease substrate.** For example, for ~20 assays, dilute 0.2 mL of EnzChek® peptidase/protease substrate (Component A) with 0.8 mL of 1X digestion buffer (from Step 1) in a disposable plastic container and mix well. Do not use glass containers. Store the solution at room temperature, protected from light, until ready to use. Table 1 is provided to facilitate the dilution of kit components for a different number of assays; alternatively, the following equations may be used:

A. Buffer Dilution. $\text{volume of 20X buffer concentrate required (mL)} = \# \text{ assays}/50$
 $\text{volume of deionized H}_2\text{O required (mL)} = (\text{volume of 20X concentrate}) \times 19$

B. Substrate Dilution. $\text{volume of 5X substrate concentrate required (mL)} = \# \text{ assays}/100$
 $\text{volume of 1X buffer required (mL)} = (\text{volume of 5X substrate concentrate}) \times 4$

4. Prepare enzyme standard and sample dilutions. Titrate concentrations of enzyme and one buffer-only control. We recommend using at least 0.5 mL of 1X digestion buffer. Detection limits for a wide variety of proteases are given in Table 2.

5. Load microplate wells. Add 50 μL of protease dilutions (from Step 4) into separate wells of a microplate, then add 50 μL of EnzChek[®] peptidase/protease substrate working solution (from Step 3) and mix well. Duplicates or triplicates are recommended. Incubate for 60 minutes at room temperature, protected from light (incubation times may vary; see “Protocol Details”).

6. Measure the fluorescence using a microplate reader. Excitation and emission maxima are 502 and 528 nm, respectively (Figure 1). Excitation/emission settings of 490/520 nm work well in the assay.

7. Use a standard curve to determine protease activity. For the protease standards, plot protease amount vs. fluorescence and fit a line to the data points. Example standard curves are given in Figures 2 and 3.

Protocol Details

Reagent Dilution

Buffers other than that provided may be required for successful use of the EnzChek[®] peptidase/protease assay. The digestion buffer in this kit is recommended for detecting the protease activity of most proteolytic enzymes with activity optima from pH 7.4 to 8.0. However, if you are working with an enzyme that requires activation compounds or a unique pH environment, prepare the specific buffer required and substitute it for the supplied digestion buffer.

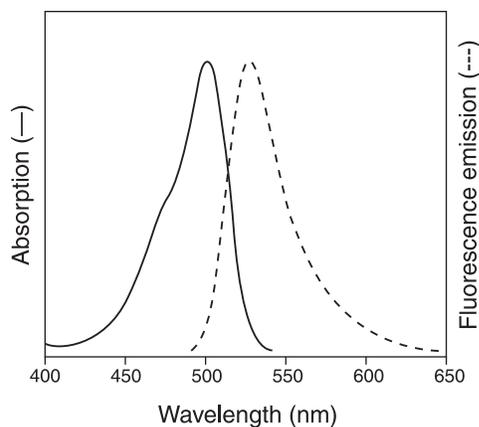


Figure 1. Normalized excitation and emission maxima for the EnzChek[®] peptidase/protease substrate digestion product in 10 mM Tris-HCl (pH 7.8).

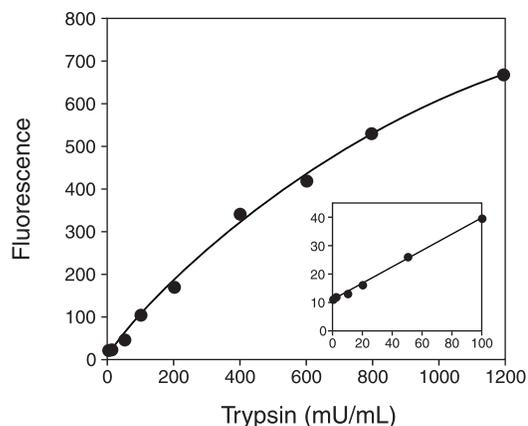


Figure 2. Sample standard curves obtained with the EnzChek[®] Peptidase/Protease Assay Kit. Trypsin (Sigma T8802, EC 3.4.21.4) was assayed in 10 mM Tris-HCl (pH 7.8) digestion buffer from 10 to 1200 mU/mL using the EnzChek[®] peptidase/protease substrate. The inset shows a separate experiment using the same enzyme, but at activities from 2 to 100 mU/mL. Samples were incubated for 60 minutes at room temperature. Fluorescence was measured at 490/520 nm; background fluorescence was subtracted from the inset data.

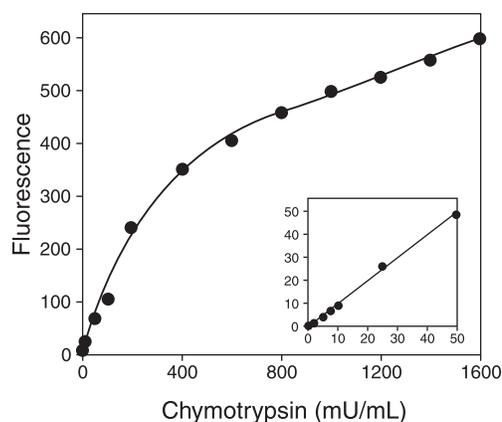


Figure 3. Sample standard curves obtained with the EnzChek[®] Peptidase/Protease Assay Kit. α -Chymotrypsin (Sigma C7762, EC 3.4.21.1) was assayed from 10 to 1600 mU/mL using the EnzChek[®] peptidase/protease substrate. The inset shows a separate experiment using the same enzyme, but at activities from 2 to 50 mU/mL. Samples were incubated for 60 minutes at room temperature. Fluorescence was measured at 490/520 nm; background fluorescence was subtracted from the inset data.

Sample Volumes

The assay has been optimized for 100 μ L total volume. We recommend preparing enzyme standard and sample dilutions in at least 0.5 mL, then aliquoting into separate microplate wells.

Assay Time and Temperature

The assay temperature is “room temperature,” defined here as 20–25°C. Assay temperatures outside of this range have not been tested, but may be acceptable. The assay described here has been optimized for a 60 minute incubation period. Sensitivity may be increased by incubating longer. Conversely, if high sensitivity is not required, incubation times may be reduced. We recommend incubations of at least 5 minutes (Figure 4). The exact time interval is not critical. However, it is important that all reactions be incubated for approximately the same time.

Standard Curves and Detection Limits

Detection limits may vary with instrumentation and enzyme source. Enzyme activity may be affected by incubation buffers and temperature, as well as the storage conditions and number of freeze–thaw cycles to which the enzyme preparation has been subjected. Different dilution schemes of the same enzyme source may also cause variation in activity. For these reasons, we recommend preparing new standard curves on the same day samples are run, preferably on the same microplate. Use an appropriate enzyme standard of known specific activity that closely matches the activity of the enzyme being assayed.

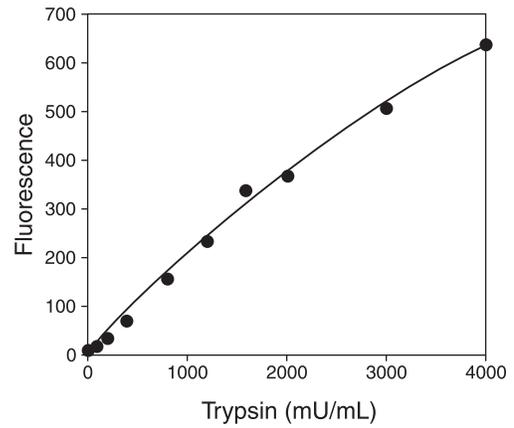


Figure 4. Incubation time with the EnzChek® Peptidase/Protease Assay Kit. Trypsin (Sigma T8802, EC 3.4.21.4) was assayed in 10 mM Tris-HCl (pH 7.8) digestion buffer from 50 to 4000 mU/mL using the EnzChek® peptidase/protease substrate. Samples were incubated for 5 minutes at room temperature and fluorescence was measured at 490/520 nm.

Table 1. Dilutions for the EnzChek® Peptidase/Protease Assay.

| Number of Assays | Buffer Dilution | | | Substrate Dilution | |
|------------------|---|---|--------------------------|-------------------------------------|-----------------------|
| | Volume of 20X Buffer (Component B) (mL) | Volume of Deionized H ₂ O (mL) | Volume 1X Available (mL) | Volume Substrate (Component A) (mL) | Volume 1X Buffer (mL) |
| 10 | 0.2 | 3.8 | 3 | 0.1 | 0.4 |
| 20 | 0.4 | 7.6 | 6 | 0.2 | 0.8 |
| 50 | 1 | 19 | 15 | 0.5 | 2 |
| 100 | 2 | 38 | 30 | 1 | 4 |

Table 2. Detection Limits of the EnzCheck® Peptidase/Protease Assay Kit.

| Protease | Source | Detection Limit (mU/mL) |
|---|-------------------------------|-------------------------|
| Trypsin from bovine pancreas | Sigma T8802 (EC 3.4.21.4) | 2 |
| α-Chymotrypsin (Type I-S) from bovine pancreas | Sigma C7762 (EC 3.4.21.1) | 2 |
| Subtilisin A from <i>Bacillus licheniformis</i> | Sigma P5380 (EC 3.4.21.62) | 0.001 |
| Protease from <i>Aspergillus oryzae</i> | Sigma P6110 (EC 3.4.21.63) | 0.004 |
| Protease from <i>Bacillus amyloliquefaciens</i> | Sigma P1236 (EC 3.4.24.28) | 1.0 × 10 ⁻⁵ |
| Protease from <i>Bacillus polymyxa</i> | Sigma P6141 (EC 3.4.24.32) | 1.5 |
| Protease from <i>Bacillus</i> sp. | Sigma P5985 (EC 3.4.21.62) | 0.005 |
| Elastase (Type IV) from porcine pancreas | Sigma E0258 (EC 3.4.21.36) | 1 |
| Thermolysin from <i>Bacillus thermoproteolyticus rokko</i> | Sigma T7902 (EC 3.4.24.27) | 1.5 |
| Papain from <i>Carica papaya</i> | Fluka 76218 (EC 3.4.22.2) | 0.08 |
| Ficin from fig tree latex | Sigma F6008 (EC 3.4.22.3) | 0.04 |
| Bromelain from pineapple stem | Sigma B4882 (EC 3.4.22.32) | 0.01 |
| Pepsin from porcine gastric mucosa | Sigma P7000 (EC 3.4.23.1) | 0.25 |

The detection limits listed here are defined as the amount of enzyme required to cause approximately a 10% change in fluorescence compared to a control sample. All samples were incubated for 60 minutes at room temperature using the EnzChek® peptidase/protease assay. Assays were performed in 10 mM Tris-HCl (pH 7.8), except for pepsin, which was assayed in 10 mM HCl (pH 1.8). Dilutions of papain, bromelain, and ficin were made from a buffer containing 30 mM L-cysteine. Enzyme unit definitions are the standard definitions for each individual enzyme. Detection limits may vary; see *Protocol Details*.

Product List Current prices may be obtained from our website or from our Customer Service Department.

| Cat # | Product Name | Unit Size |
|--------|--|-----------|
| E33758 | EnzChek® Peptidase/Protease Assay Kit *100 assays* | 1 kit |

Contact Information

Further information on Molecular Probes products, including product bibliographies, is available from your local distributor or directly from Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Paisley, United Kingdom. All others should contact our Technical Service Department in Eugene, Oregon.

Please visit our website — probes.invitrogen.com — for the most up-to-date information.

Molecular Probes, Inc.

29851 Willow Creek Road, Eugene, OR 97402
Phone: (541) 465-8300 • Fax: (541) 335-0504

Customer Service: 6:00 am to 4:30 pm (Pacific Time)

Phone: (541) 335-0338 • Fax: (541) 335-0305 • probesorder@invitrogen.com

Toll-Free Ordering for USA:

Order Phone: (800) 438-2209 • Order Fax: (800) 438-0228

Technical Service: 8:00 am to 4:00 pm (Pacific Time)

Phone: (541) 335-0353 • Toll-Free (800) 438-2209
Fax: (541) 335-0238 • probestech@invitrogen.com

Invitrogen European Headquarters

Invitrogen, Ltd.
3 Fountain Drive
Inchinnan Business Park
Paisley PA4 9RF, UK
Phone: +44 (0) 141 814 6100 • Fax: +44 (0) 141 814 6260
Email: euroinfo@invitrogen.com
Technical Services: eurotech@invitrogen.com

Molecular Probes products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

Limited Use Label License

For research use only. Not intended for any animal or human therapeutic or diagnostic use. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402. Tel: (541) 465-8300. Fax: (541) 335-0504.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation ® are registered with the U.S. Patent and Trademark Office.

Copyright 2005, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.