



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: info@bpsbioscience.com

**Data sheet**  
***Cathepsin B Inhibitor Screening Assay Kit***  
**Catalog #79590**  
**Size: 96 reactions**

**BACKGROUND:** Cathepsin B is a cysteine protease that primarily functions as an endopeptidase within endolysosomal compartments in normal cells. High levels of cathepsin B are found in a wide variety of human cancers and in experimental models, such as transgenic models of murine pancreatic and mammary carcinomas, causal roles for this protein have been demonstrated in initiation, growth/tumor cell proliferation, angiogenesis, invasion, and metastasis.

**DESCRIPTION:** The *Cathepsin B Inhibitor Screening Assay Kit* is designed to measure the protease activity of Cathepsin B for screening and profiling applications. The Cathepsin B assay kit comes in a convenient 96-well format, with purified Cathepsin B, its fluorogenic substrate, and Cathepsin buffer for 100 enzyme reactions. In addition, the kit includes the cathepsin inhibitor E-64 for use as a control inhibitor.

**COMPONENTS:**

Catalog #	Component	Amount	Storage	
80001	Cathepsin B	10 µg	-80 °C	Avoid multiple freeze/thaw cycles!
80349	Fluorogenic Cathepsin Substrate 1 (5 mM)	10 µl	-20 °C	
	4X Cathepsin buffer	2 ml	-20 °C	
	E-64 (1 mM)	10 µl	-20 °C	
	96-well black microplate			

**MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

0.5 M DTT in aqueous solution  
Adjustable micropipettor and sterile tips  
Fluorescent microplate reader

**APPLICATIONS:** Great for studying enzyme kinetics and screening small molecular inhibitors for drug discovery and HTS applications.

**STABILITY:** Up to 6 months from date of receipt, when stored as recommended.

**REFERENCES:**

1. Aggarwal, N., & Sloane, B. F. (2014). Cathepsin B: multiple roles in cancer. *PROTEOMICS–Clinical Applications* **8(5-6)**, 427-437.
2. Siklos, M., BenAissa, M., & Thatcher, G. R. (2015). Cysteine proteases as therapeutic targets: does selectivity matter? A systematic review of calpain and cathepsin inhibitors. *Acta Pharmaceutica Sinica B*, **5(6)**, 506-519.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: info@bpsbioscience.com

#### ASSAY PROTOCOL:

*All samples and controls should be tested in duplicate.*

- 1) Prepare **1X Cathepsin buffer** by diluting **4X Cathepsin buffer** 4-fold and 0.5 M DTT 100-fold (5 mM final assay concentration) into water. For example, to prepare 10 ml, add 2.5 ml of **4X Cathepsin buffer** and 0.1 ml of 0.5 M DTT to 7.4 ml of water.
- 2) Add 20  $\mu$ l **1X Cathepsin buffer** to each well designated "Blank".
- 3) Thaw **Cathepsin B** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Activate **Cathepsin B** by diluting the enzyme to 10 ng/ $\mu$ l in **1X Cathepsin buffer** for 15 minutes at room temperature. Aliquot remaining **Cathepsin B** enzyme into single-use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. Note: **Cathepsin B** is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
- 4) Prepare **Enzyme solution** (0.02 ng/ $\mu$ l Cathepsin B) by diluting **10 ng/ $\mu$ l diluted, activated Cathepsin B** 500-fold in **1X Cathepsin buffer**. For example, to prepare 1000  $\mu$ l, add 2  $\mu$ l **10 ng/ $\mu$ l activated Cathepsin B** to 998  $\mu$ l **1X Cathepsin buffer**.
- 5) Add 20  $\mu$ l **Enzyme solution** (0.02 ng/ $\mu$ l Cathepsin B) to each well designated "Positive Control", "Negative Control" and "Test Inhibitor".
- 6) Add 5  $\mu$ l **Inhibitor solution** to each well designated "Test Inhibitor". For the wells labeled "Positive Control" and "Blank", add 5  $\mu$ l **Inhibitor buffer** (same solution without inhibitor). Optional: for each well designated "Negative Control", add 5  $\mu$ l **E-64** diluted 0.1 – 0.0001  $\mu$ M in **1X Cathepsin buffer**. Incubate at room temperature for 10 minutes.
- 7) Prepare **Substrate solution** (10  $\mu$ M) by diluting **Fluorogenic Cathepsin Substrate 1** (5 mM) 500-fold in **1X Cathepsin buffer**. For example, to prepare 1000  $\mu$ l, add 2  $\mu$ l **Fluorogenic Cathepsin Substrate 1** (5 mM) to 998  $\mu$ l **1X Cathepsin buffer**. Store remaining undiluted substrate in aliquots at -20°C. Do not re-use diluted substrate.
- 8) Add **25  $\mu$ l Substrate solution** (10  $\mu$ M) to all wells. Incubate reaction at room temperature for 60 minutes.

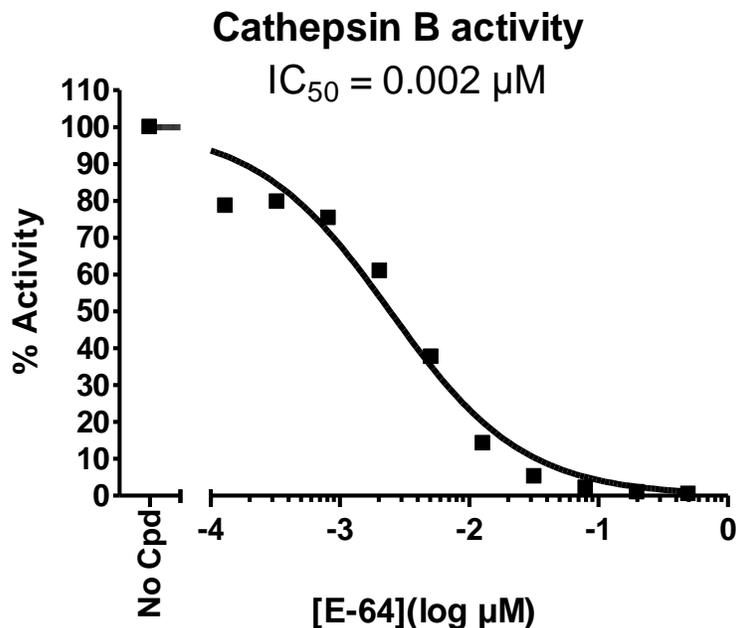
	Positive Control	Negative Control	Test Inhibitor	Blank
Enzyme solution (0.02 ng/ $\mu$ l Cathepsin B)	20 $\mu$ l	20 $\mu$ l	20 $\mu$ l	-
1X Cathepsin buffer	-	-	-	20 $\mu$ l
Inhibitor (in Cathepsin buffer)	-	-	5 $\mu$ l	-
Inhibitor buffer (no inhibitor)	5 $\mu$ l	-	-	5 $\mu$ l
E-64	-	5 $\mu$ l	-	
Substrate solution (10 $\mu$ M)	25 $\mu$ l	25 $\mu$ l	25 $\mu$ l	25 $\mu$ l
Total	50 $\mu$ l	50 $\mu$ l	50 $\mu$ l	50 $\mu$ l

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

- 9) Read fluorescence intensity of the samples ( $\lambda_{\text{excitation}} = 360 \text{ nm}$ ;  $\lambda_{\text{emission}} = 460 \text{ nm}$ ) in an appropriate microplate reader. "Blank" value is subtracted from all readings.

**Example of assay results:**



Cathepsin B inhibition by E-64, measured using the *Cathepsin B Inhibitor Screening Assay Kit*, BPS Bioscience, #79590. Fluorescence was measured using a Bio-Tek microplate reader. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**RELATED PRODUCTS:**

<u>Product</u>	<u>Catalog#</u>	<u>Size</u>
Cathepsin B	80001	10 $\mu\text{g}$
Cathepsin L	80005	10 $\mu\text{g}$
Cathepsin S	80008	10 $\mu\text{g}$
Cathepsin V	80009	10 $\mu\text{g}$
Fluorogenic Cathepsin Substrate 1	80349	100 $\mu\text{l}$
Fluorogenic Cathepsin F Substrate	80350	100 $\mu\text{l}$

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)