

Description

The KLK11 Fluorogenic Assay Kit is designed to measure KLK11 (Kallikrein-11) protease activity for screening and profiling applications. The assay kit comes in a convenient 96-well format, with enough recombinant KLK11 (amino acids 54-282), fluorogenic AMC-based peptide substrate, and assay buffer for 100 enzyme reactions.

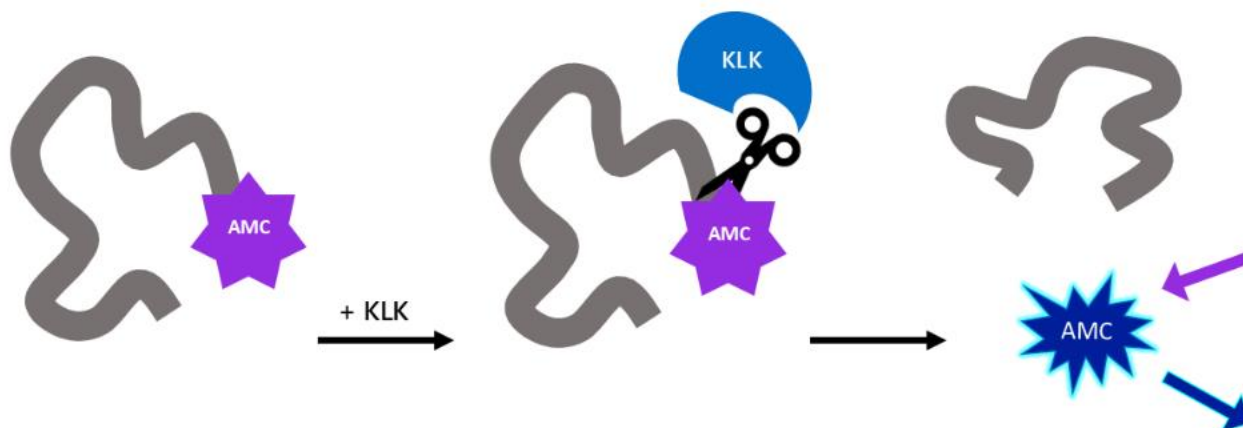


Figure 1: Illustration of the assay principle.

A peptide substrate conjugated to AMC is incubated with KLK11. The conjugated AMC displays only weak fluorescence. Upon proteolysis, AMC is no longer quenched and exhibits strong blue fluorescence. The increase in fluorescence is proportional to the KLK11 activity.

Background

KLK11 (kallikrein 11), also known as TLSP (Trypsin-like serine protease), is a secreted serine endopeptidase that participates in skin desquamation. Abnormal expression of this protein can result in skin barrier related disorders. It can be found in multiple tissues, such as the brain, skin, breast, prostate and intestine. It can be found at high levels in the serum of breast and prostate cancer patients, and it is considered a favorable prognostic factor. However, in other cancer types it seems to correlate with a poor prognosis. In a model of colorectal cancer (CRC) the knockdown of KLK11 lead to inhibition of cell growth and increased in apoptosis. A deeper understanding of the role in health and disease of this protein may open new therapeutic opportunities in oncology.

Applications

Enzyme kinetics studies and screening small molecule inhibitors for drug discovery and high-throughput screening (HTS) applications.

Supplied Materials

Catalog #	Name	Amount	Storage
84280-KC12 ⁺	Kallikrein-11(KLK11), Lyophilized	12 µg	-80°C
84003-KC5	PR Substrate 4	5 µl	-80°C
84004-KC10	PR-03 Assay Buffer	10 ml	-80°C
79685	96-well black microplate	1	Room Temp

Materials Required but Not Supplied

- Adjustable micropipettor and sterile tips
- Fluorimeter capable of excitation at $\lambda=380$ nm and detection at $\lambda=460$ nm

Stability



This assay kit will perform optimally for up to **6 months** from date of receipt when the materials are stored as directed.

Safety



This product is for research purposes only and not for human or therapeutic use. This product should be considered hazardous and is harmful by inhalation, in contact with skin, eyes, clothing, and if swallowed. If contact occurs, wash thoroughly.

Contraindications

- The final concentration of DMSO in the assay should not exceed 1%.
- Compounds that are fluorescent may interfere with the results, depending on their spectral excitation and emission properties.
- It is recommended that the compound alone is tested to determine any potential interference of the compound with the assay results.

Assay Protocol

- All samples and controls should be performed in duplicates.
 - The assay should include “Negative Control”, “Positive Control,” and a “Test Inhibitor” conditions.
 - We recommend maintaining the diluted protein on ice during use.
 - For detailed information on protein handling please refer to [Protein FAQs \(bpsbioscience.com\)](https://bpsbioscience.com/protein-faqs/).
 - We recommend using Gabexate Mesylate (BPS Bioscience #84177) as internal control. If not running a dose response curve for the control inhibitor, we recommend running the control inhibitor at 0.1X, 1X and 10X the IC₅₀ value shown in the validation data below.
1. Resuspend KLK11 with 24 μ l of distilled water to get a concentration of 0.5 mg/ml.
 2. Dilute 100-fold the **KLK11** to 5 ng/ μ l with **PR-03 Assay Buffer** (you need 20 μ l/well).
 3. Add 20 μ l of **diluted KLK11** to all wells except “Negative Control” wells.
 4. Add 20 μ l of **PR-03 Assay Buffer** to the “Negative Control” wells.
 5. Prepare the **Test inhibitor** (5 μ l/well): for a titration, prepare serial dilutions at concentrations 10-fold higher than the desired final concentrations. The final volume of the reaction is 50 μ l.
 - 5.1. If the Test Inhibitor is water-soluble, prepare serial dilutions in the PR-03 Assay Buffer at concentrations 10-fold higher than the desired final concentrations.

For positive and negative controls, use PR-03 Assay Buffer (Diluent Solution).

OR

5.2. If the Test inhibitor is soluble in DMSO, prepare the test inhibitor in 100% DMSO at a concentration 100-fold higher than the highest desired final concentration, then dilute the inhibitor 10-fold in PR-03 Assay Buffer to prepare the highest concentration of the serial dilutions. The concentration of DMSO is now 10%.

Using PR-03 Assay Buffer containing 10% DMSO to keep the concentration of DMSO constant, prepare serial dilutions of the Test Inhibitor at 10-fold the desired final concentrations.

For positive and negative controls, prepare 10% DMSO in PR-03 Assay Buffer (vol/vol) so that all wells contain the same amount of DMSO (Diluent Solution).

Note: The final concentration of DMSO in the assay should not exceed 1%.

6. Add 5 µl of **Test inhibitor** to each well designated “Test Inhibitor”.
7. Add 5 µl of **Diluent Solution** to the “Positive Control” and “Negative Control” wells.
8. Preincubate the Test inhibitor with the diluted KLK11 for 30 minutes at Room Temperature (RT) with gentle agitation.
9. Dilute 500-fold the **PR Substrate 4** with PR-03 Assay Buffer (25 µl/ well).
10. Initiate the reaction by adding 25 µl of **diluted PR Substrate 4** to all wells.



Protect your samples from direct exposure to light.

Component	Negative Control	Positive Control	Test Inhibitor
PR-03 Assay Buffer	20 µl	-	-
Test Inhibitor	-	-	5 µl
Diluent Solution	5 µl	5 µl	-
Diluted KLK11 (5 ng/µl)	-	20 µl	20 µl
Pre-incubate 30 minutes at Room Temperature			
Diluted PR Substrate 4	25 µl	25 µl	25 µl
Total	50 µl	50 µl	50 µl

11. Incubate at RT for 2 hours.
12. Read the fluorescence intensity of the samples ($\lambda_{\text{excitation}} = 380 \text{ nm}$; $\lambda_{\text{emission}} = 460 \text{ nm}$, bandwidth = 10) in a fluorescence reader.

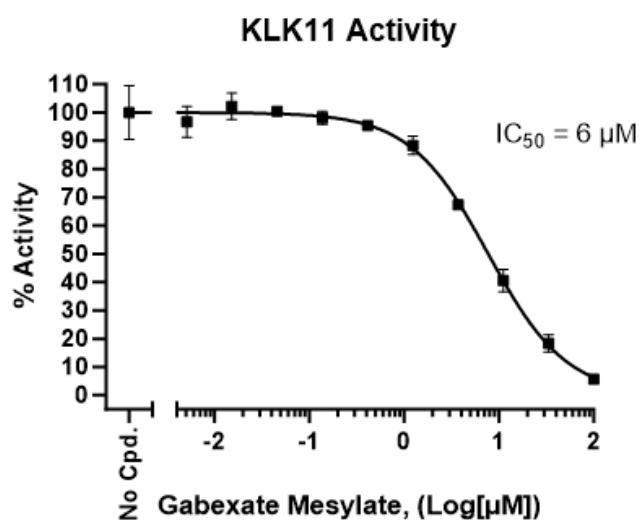
Example Results:

Figure 2. KLK11 activity is inhibited by Gabexate Mesylate.

KLK11 activity was measured in the presence of increasing concentrations of Gabexate Mesylate (#84177). Results are expressed as percentage of activity relative to the positive control (measured in the absence of inhibitor and set at 100%).

Data shown is representative.

Troubleshooting Guide

Visit bpsbioscience.com/assay-kits-faq for detailed troubleshooting instructions. For lot-specific information and all other questions, please visit <https://bpsbioscience.com/contact>.

References

- Li Y.-C., *et al.*, 2022 *Biomedicines* 10 (2): 474.
 Xu Z., *et al.*, 2016 *Exp Ther Med* 12(5): 2855-2860.

Version 062526