

# TACE Inhibitor Assay Kit ab155889

[2 Images](#)

### Overview

|                         |  |
|-------------------------|--|
| <b>Product name</b>     | TACE Inhibitor Assay Kit   |
| <b>Detection method</b> | Fluorescent  |
| <b>Assay type</b>       | Quantitative   |
| <b>Assay time</b>       | 0h 20m   |
| <b>Product overview</b> | In Abcam's TACE inhibitor screening Kit, TACE (tumor necrosis factor- $\alpha$ -converting enzyme) hydrolyzes the specific FRET substrate to release the quenched fluorescent group, which can be detected fluorometrically at Ex/Em = 318/449 nm. In the presence of the potent TACE inhibitor, the hydrolyzation of substrate will be impeded. The kit provides a rapid, simple, sensitive and reliable test suitable for high-throughput screening of TACE inhibitors and can be modified to check the relative TACE activity. Inhibitor Control GM6001 is included to compare the efficacy of test inhibitors. |

|              |   |
|--------------|---|
| <b>Notes</b> | <p>This product is manufactured by BioVision, an Abcam company and was previously called K366 TACE Inhibitor Screening Assay Kit ( Fluorometric). K366-100 is the same size as the 100 test size of ab155889.</p> <p>The TACE (tumor necrosis factor-<math>\alpha</math>-converting enzyme), also called ADAM metallopeptidase domain 17 (ADAM17), is a 70-kDa enzyme that belongs to the ADAM protein family of disintegrins and metalloproteases. TACE is believed to be involved in the processing of tumor necrosis factor alpha (TNF-<math>\alpha</math>) at the surface of the cell, and from within the intracellular membranes of the trans-Golgi network. This process, which is also known as 'shedding', involves the cleavage and release of a soluble ectodomain from membrane-bound pro-proteins (such as pro-TNF-<math>\alpha</math>), and is of known physiological importance.</p> <p>The standard is affinity purified.</p> |
|--------------|---|

|                 |                   |
|-----------------|-------------------|
| <b>Platform</b> | Microplate reader |
|-----------------|-------------------|

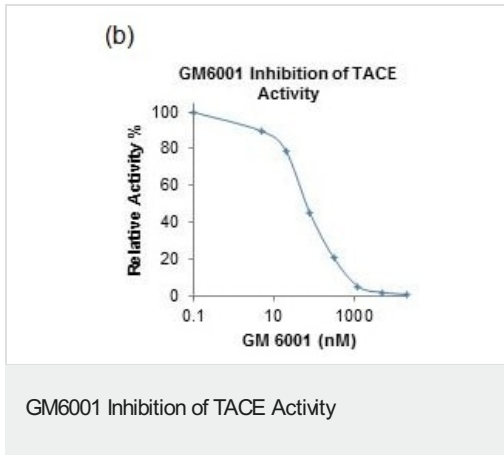
### Properties

|                             |  |
|-----------------------------|--|
| <b>Storage instructions</b> | Store at -20°C. Please refer to protocols. |
|-----------------------------|--|

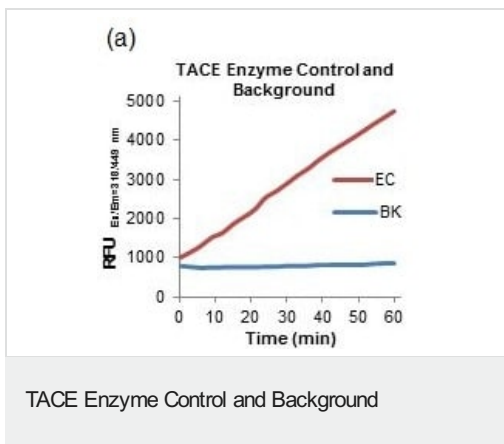
| Components        | 100 tests |
|-------------------|-----------|
| TACE Assay Buffer | 1 x 25ml  |
|                   |           |

| Components             | 100 tests |
|------------------------|-----------|
| TACE Enzyme            | 1 vial    |
| TACE Inhibitor Control | 1 x 20µl  |
| TACE Substrate         | 1 x 0.2ml |

## Images



TACE activity with increasing GM6001 inhibition. Produced following kit protocol.



Fluorescence of TACE enzyme control and background. Produced following kit protocol.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors