



6405 Mira Mesa Blvd Ste 100  
San Diego Ca 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

**Data Sheet**  
**ADAM17 Fluorogenic Assay Kit**  
Catalog #78000  
Size: 96 reactions

**BACKGROUND:** ADAM17 (ADAM Metallopeptidase Domain 17) is part of the ADAM family of disintegrins and metalloproteases. Initially identified as TNF- $\alpha$  converting enzyme (TACE), ADAM 17 has been linked to a number of diverse signaling pathways. It cleaves ectodomains of various transmembrane proteins and regulates cytokine shedding. It also plays a role in inflammatory skin and bowel disease.

**DESCRIPTION:** The *ADAM17 Fluorogenic Assay Kit* is provided in a convenient 96-well format, with purified ADAM17, ADAM Fluorogenic Substrate, and ADAM assay buffer for 96 enzyme reactions. The key to the *ADAM17 Fluorogenic Assay Kit* is the fluorogenic substrate. Using this kit, only one simple step on a microtiter plate is required for ADAM17 reactions. A sample containing ADAM17 is incubated in a reaction mixture with the fluorogenic substrate and fluorescence ( $\lambda_{ex} = 485 \text{ nm}$ ,  $\lambda_{em} = 530 \text{ nm}$ ) is measured using a plate reader.

**COMPONENTS:**

Catalog #	Component	Amount	Storage	
	ADAM17	3 $\mu\text{g}$	-80°C	<b>Avoid freeze/thaw cycles!</b>
	ADAM Fluorogenic Substrate (1 mM)	50 $\mu\text{l}$	-80°C	
78001	1X ADAM Assay Buffer	5 ml	-20°C	
79685	96-well black plate	1	Room temp.	

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

**STABILITY:** One year from date of receipt when stored as directed.

**REFERENCES:**

- Scheller, Jürgen, *et al.* 2011. "ADAM17: a molecular switch to control inflammation and tissue regeneration." *Trends in Immunology* **32(8)**: 380-387.
- Blaydon, Diana C., *et al.* 2011. "Inflammatory skin and bowel disease linked to ADAM17 deletion." *New England Journal of Medicine* **365(16)**: 1502-1508.

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

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**  
Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd Ste 100  
San Diego Ca 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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

Luminometer or Fluorogenic microplate reader capable of reading chemiluminescence  
Adjustable micropipettor and sterile tips  
Rotating or rocker platform

**ASSAY PROTOCOL:**

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

**Step 1:**

- 1) Prepare the master mixture: N wells x (24.5  $\mu$ l **1x ADAM Assay Buffer 1** + 0.5  $\mu$ l **ADAM Fluorogenic Substrate** (1 mM)).
- 2) Add 25  $\mu$ l of master mixture to each well designated for the "Positive Control," "Test Inhibitor," and "Blank."
- 3) Thaw **ADAM17** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **ADAM17** into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. *Note: ADAM17 enzyme is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 4) Dilute **ADAM17** in **1X ADAM Assay Buffer** at 1.25 ng/ $\mu$ l (25 ng/reaction). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.
- 5) Prepare the test inhibitor solution.

The final concentration of DMSO in the assay should not exceed 1%. If the inhibitor compound is dissolved in DMSO, make a 100-fold higher concentration of the compound than the highest concentration you want to test in DMSO. Then make a 10-fold dilution in **1X ADAM Assay Buffer** (at this step the compound concentration is 10-fold higher than the final concentration in 10% DMSO). To determine an IC<sub>50</sub> or to test lower concentrations of the compound, prepare a series of further dilutions in **1X ADAM Assay Buffer** containing 10% DMSO (the final concentration of the DMSO will be 1% in all samples).

If the inhibitor compound is dissolved in water, make a solution of the compound 10-fold higher than the final concentration in **1X ADAM Assay Buffer**.

- 6) Add 5  $\mu$ l of test Inhibitor solution to each well labeled as "Test Inhibitor." For the "Positive Control" and "Blank," add 5  $\mu$ l of inhibitor buffer (same solution without inhibitor compound; usually 10% DMSO in **1X ADAM Assay Buffer**).
- 7) Add 20  $\mu$ l of **1X ADAM Assay Buffer** to the wells designated "Blank."

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

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

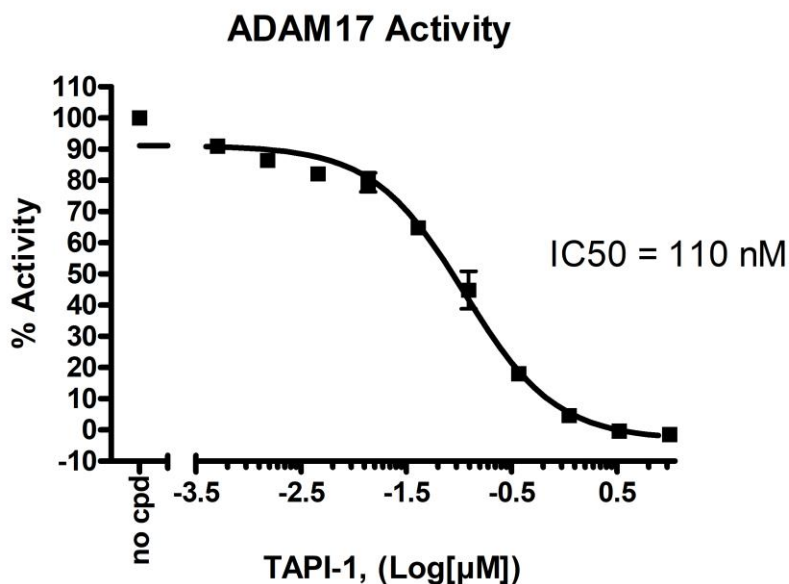
- 8) Initiate reaction by adding 20  $\mu\text{l}$  of diluted **ADAM17** (1.25 ng/ $\mu\text{l}$ ) to the wells designated "Positive Control" and "Test Inhibitor." Incubate for 60 minutes at room temperature.

	Positive Control	Test Inhibitor	Blank
Master Mixture	25 $\mu\text{l}$	25 $\mu\text{l}$	25 $\mu\text{l}$
Test Inhibitor	-	5 $\mu\text{l}$	-
Inhibitor Buffer	5 $\mu\text{l}$	-	5 $\mu\text{l}$
1x ADAM Assay Buffer	-	-	20 $\mu\text{l}$
ADAM17 (1.25 ng/ $\mu\text{l}$ )	20 $\mu\text{l}$	20 $\mu\text{l}$	-
<b>Total</b>	<b>50 <math>\mu\text{l}</math></b>	<b>50 <math>\mu\text{l}</math></b>	<b>50 <math>\mu\text{l}</math></b>

### Step 2:

- 1) Read fluorescence at  $\lambda_{\text{ex}} = 485 \text{ nm}$ ,  $\lambda_{\text{em}} = 530 \text{ nm}$ . "Blank" value is subtracted from all measurements.

### Example of Assay Results:



ADAM17 enzyme activity, measured using the *ADAM17 Assay Kit*, BPS Bioscience #78000. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd Ste 100  
 San Diego Ca 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

### TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Fluorogenic signal of positive control reaction is weak	ADAM17 enzyme has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh enzyme. Store enzyme in single-use aliquots. Increase time of enzyme incubation. Increase enzyme concentration.
	Incorrect wavelength	Use correct filters for Fluorogenic plate reader. Reading at wavelengths outside of 455 ± 10 nm will give a decreased signal.
Fluorogenic signal is erratic or varies widely among wells	Inaccurate pipetting/technique	Run duplicates of all reactions. Use a multichannel pipettor. Use master mixes to minimize errors.
	Bubbles in wells	Pipette slowly to avoid bubble formation. Tap plate lightly to disperse bubbles; be careful not to splash between wells.
Background (signal to noise ratio) is high	Sample solvent is inhibiting the enzyme	Run negative control assay including solvent. Maintain DMSO level at <1% Increase time of enzyme incubation.
	Results are outside the linear range of the assay	Use different concentrations of ADAM17 enzyme to create a standard curve.

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

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**  
 Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)