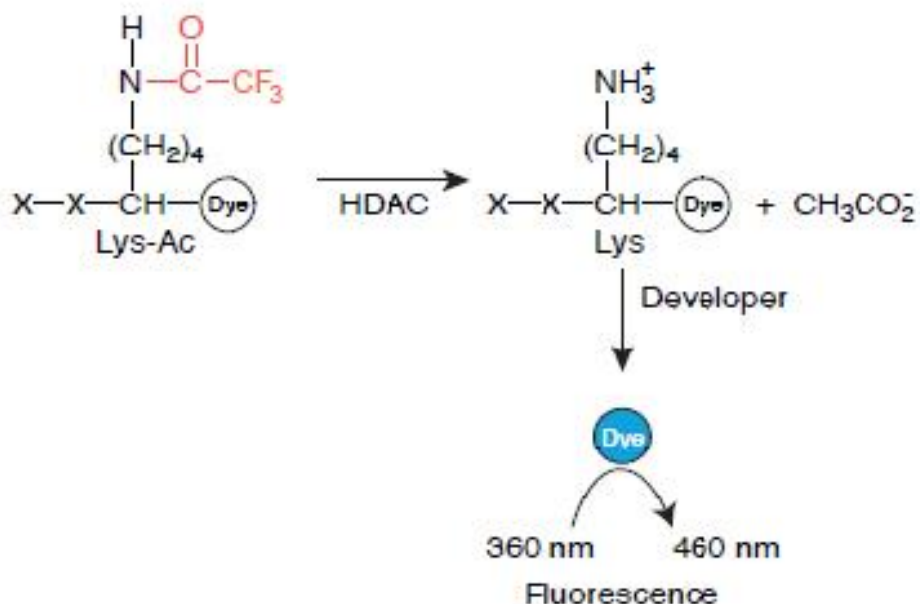


## Data Sheet

### **HDAC11 Fluorogenic Assay Kit** Catalog #: 50687

**DESCRIPTION:** The *HDAC11 Fluorogenic Assay Kit* is a complete assay system designed to measure histone deacetylase 11 (HDAC11) activity for screening and profiling applications. It comes in a convenient 96-well format, with all the reagents necessary for 100 fluorescent HDAC11 activity measurements. In addition, the kit includes purified HDAC11 enzyme and a potent HDAC inhibitor, Trichostatin A, for use as a positive and negative control. The *HDAC11 Fluorogenic Assay Kit* is based on a unique fluorogenic substrate and developer combination. This assay method eliminates dealing with the radioactivity, extraction, and chromatography aspects of traditional assays. Using this kit, only two simple steps on a microtiter plate are needed to analyze the HDAC11 activity level. First, the HDAC fluorometric substrate, containing an acetylated lysine side chain, is incubated with purified HDAC11. The deacetylation sensitizes the substrate so subsequent treatment with the Lysine Developer produces a fluorophore that can then be measured using a fluorescence reader.

HDACs regulate cellular processes by catalyzing the hydrolysis of an acetyl group from acetyllysines in modified proteins. In the HDAC assay, fluorescent-dye molecules are attached to a peptide containing acetyllysine. Attachment to the peptide quenches the fluorescence of the dye. After treatment of the peptide with an HDAC, the reaction is mixed with a development solution that is specific for nonacetylated lysines. If the acetyl group has been removed from the lysine by the HDAC, this solution will release the dye allowing for fluorescence. The amount of fluorescence is therefore directly related to the level of HDAC activity.



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: [info@bpsbioscience.com](mailto:info@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

#### COMPONENTS:

| Cat. # | Component   | Amount  | Storage    |   |
|--------|---|---------|------------|---|
| 50021  | HDAC11 human recombinant enzyme                       | 10 µg   | -80°C      | <b>Avoid<br/>freeze/<br/>thaw<br/>cycles!</b> |
| 50040  | Fluorogenic HDAC substrate class 2A (5 mM)            | 25 µl   | -80°C      |   |
| 50030  | 2x HDAC Developer (contains Trichostatin A)<br>(2 µM) | 6 ml    | -80°C      |   |
|        | Trichostatin A (1 mM) in DMSO                         | 100 µl  | -20°C      |   |
| 79320  | HDAC11 Assay Buffer                                   | 10 ml   | -20°C      |   |
| 79685  | black, low binding NUNC black microtiter plate        | 1 plate | Room temp. |   |

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

0.1% solution (1 mg/ml) of bovine serum albumin (BSA) in water

Fluorimeter capable of excitation at 350-380 nm and detection at 440-460 nm

Adjustable micropipettor and sterile tips

Rotating or rocker platform

**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.

**REFERENCE:** Deubzer, H.E., *et al.*, *Int. J. Cancer*. 2013 May 1; **132(9)**:2200-8.

#### ASSAY PROTOCOL:

##### Immediately prior to assay:

- 1) Dilute **Trichostatin A** 1 mM stock 10-fold with **HDAC11 Assay Buffer** to make a 100 µM solution. Make only sufficient quantity needed for the assay; store remaining 1 mM **Trichostatin A** stock solution in aliquots at -80°C.
- 2) Dilute **HDAC substrate** 5 mM stock 250-fold with **HDAC11 Assay Buffer** to make a 20 µM solution. Make only sufficient quantity needed for the assay; store remaining 5 mM stock solution in aliquots at -80°C.
- 3) Dilute **HDAC11** in **HDAC11 Assay Buffer** to 20 ng/µl (100 ng/reaction)\*. Aliquot any remaining enzyme and store undiluted at -80°C. Keep diluted enzyme on ice. Discard any remaining diluted enzyme after use. *\*Note: optimal enzyme concentration may vary with the specific activity of the enzyme.*

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: [info@bpsbioscience.com](mailto:info@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

### Step 1:

In duplicate, add the reaction mixtures (below) to the microtiter black plate as follows:

- 1) Prepare the master mixture: N wells × (5 µl **HDAC substrate** (20 µM) + 5 µl BSA (1 mg/ml) + 30 µl **HDAC11 Assay Buffer**). Add 40 µl of master mixture to all wells.
- 2) Add 5 µl of inhibitor solution of each well designated "Test Inhibitor." For the "Positive Control" and "Blank," add 5 µl of the same solution without inhibitor (inhibitor buffer). Add 5 µl of diluted **Trichostatin A** (100 µM) to the wells designated "Inhibitor Control." Keep final DMSO concentration at or below 1%.
- 3) Add 5 µl of **HDAC11 Assay Buffer** to the wells designated "Blank."
- 4) Initiate reaction by adding 5 µl of diluted **HDAC11 enzyme** to the wells designated "Positive Control," "Test Inhibitor," and "Inhibitor Control." Incubate at 37°C for 30 minutes.

|                                 | Blank        | Positive Control | Test Inhibitor | Inhibitor Control |
|---------------------------------|--------------|------------------|----------------|-------------------|
| HDAC substrate (20 µM)          | 5 µl         | 5 µl             | 5 µl           | 5 µl              |
| BSA (1 mg/ml)                   | 5 µl         | 5 µl             | 5 µl           | 5 µl              |
| HDAC11 Assay Buffer             | 35 µl        | 30 µl            | 30 µl          | 30 µl             |
| Diluted Trichostatin A (100 µM) | -            | -                | -              | 5 µl              |
| Test Inhibitor                  | -            | -                | 5 µl           | -                 |
| Inhibitor buffer (no inhibitor) | 5 µl         | 5 µl             | -              | -                 |
| Diluted HDAC11 (20 ng/µl)       | -            | 5 µl             | 5 µl           | 5 µl              |
| <b>Total</b>                    | <b>50 µl</b> | <b>50 µl</b>     | <b>50 µl</b>   | <b>50 µl</b>      |

### Step 2:

Add 50 µl of undiluted **HDAC Developer (2x)** to each well. Incubate the plate at room temperature for 15 minutes.

### Step 3:

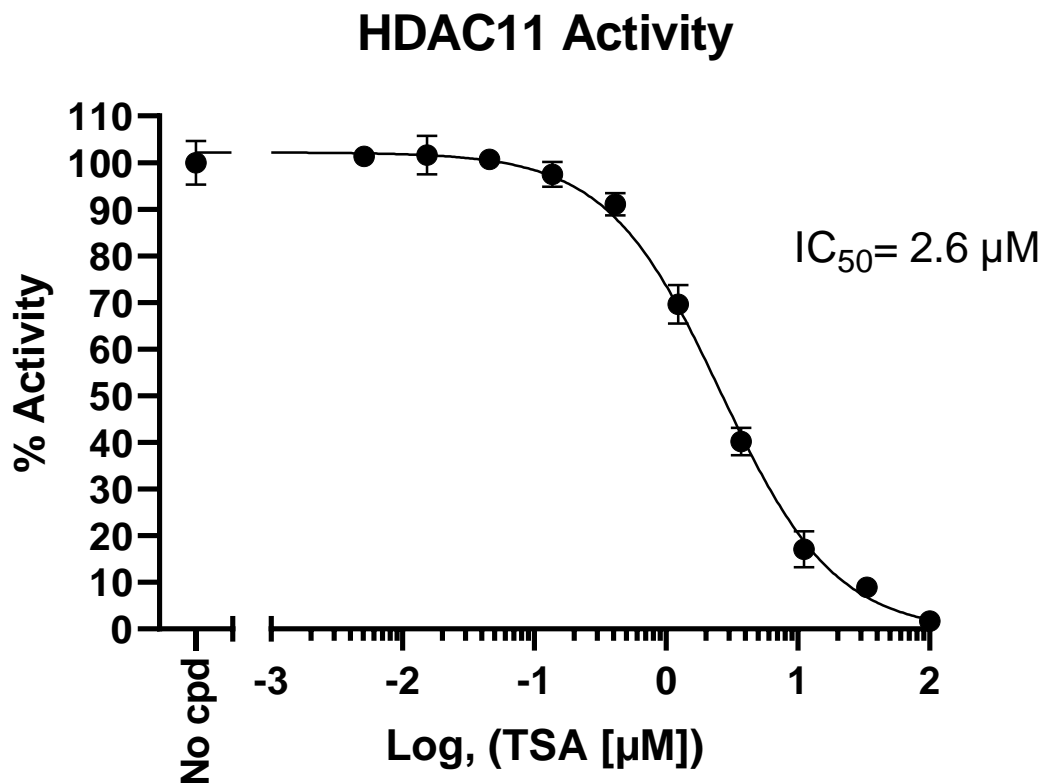
Read sample in a microtiter plate-reading fluorimeter capable of excitation at a wavelength in the range of 350-380 nm and detection of emitted light in the range of 440-460 nm. "Blank" value is subtracted from all other values.

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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Example of Assay Results:



Inhibition of HDAC11 enzyme activity by TSA, measured using the *Fluorogenic HDAC11 Assay Kit*, BPS Bioscience Catalog #50687. Fluorescence was measured using a Tecan Infinite M1000 microplate reader. *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: [info@bpsbioscience.com](mailto:info@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)

**RELATED PRODUCTS:**

| <b><u>Product Name</u></b> | <b><u>Catalog #</u></b> | <b><u>Size</u></b> |
|----------------------------|-------------------------|--------------------|
| HDAC1                      | 50051                   | 50 µg              |
| HDAC2 (C-His)              | 50002                   | 50 µg              |
| HDAC2 (C-Flag)             | 50052                   | 50 µg              |
| HDAC3/NcoR2                | 50003                   | 50 µg              |
| HDAC11                     | 50004                   | 10 µg              |
| HDAC5                      | 50005                   | 10 µg              |
| HDAC6 (C-Flag)             | 50056                   | 50 µg              |
| HDAC6 (N-GST)              | 50006                   | 50 µg              |
| HDAC6 (H216A)              | 50046                   | 50 µg              |
| HDAC6 (H611A)              | 50066                   | 50 µg              |
| HDAC7                      | 50007                   | 10 µg              |
| HDAC8                      | 50008                   | 50 µg              |
| HDAC9                      | 50009                   | 10 µg              |
| HDAC10                     | 50010                   | 50 µg              |
| HDAC11                     | 50011                   | 50 µg              |
| HDAC Assay Kit             | 50033                   | 96 reactions       |
| HDAC Assay Kit (Green)     | 50034                   | 96 reactions       |
| HDAC Class 2a Assay Kit    | 50041                   | 96 reactions       |
| HDAC2 Assay Kit            | 50062                   | 96 reactions       |
| HDAC3 Assay Kit            | 50073                   | 96 reactions       |
| HDAC5 Assay Kit            | 50065                   | 96 reactions       |
| HDAC6 Assay Kit            | 50076                   | 96 reactions       |
| HDAC7 Assay Kit            | 50077                   | 96 reactions       |
| HDAC8 Assay Kit            | 50068                   | 96 reactions       |
| HDAC9 Assay Kit            | 50069                   | 96 reactions       |
| Trichostatin A             | 27102                   | 5 mg               |
| SAHA (Vorinostat)          | 27006                   | 100 mg             |
| 4-iodo-SAHA                | 27205                   | 250 nmol           |
| Tubastatin A HCl           | 27108                   | 10 mg              |

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: [\*\*info@bpsbioscience.com\*\*](mailto:info@bpsbioscience.com)