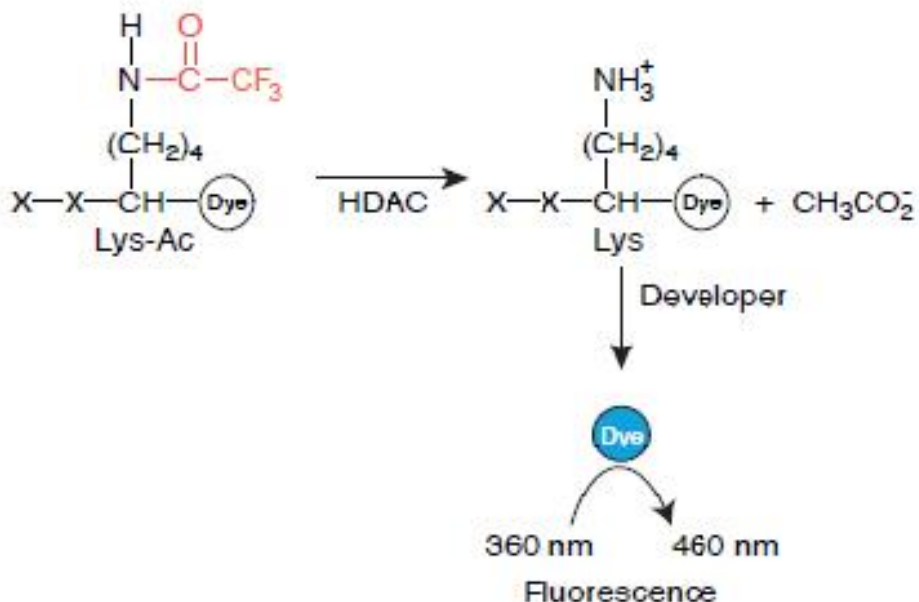


Data Sheet

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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

COMPONENTS:

Cat. #	Component	Amount	Storage	
50021	HDAC11 human recombinant enzyme	10 µg	-80°C	Avoid freeze/ thaw cycles!
50040	Fluorogenic HDAC substrate class 2A (5 mM)	50 µl	-80°C	
50030	2x HDAC Developer (contains Trichostatin A) (50 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 10 ng/µl (50 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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@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 \times (5 μ l **HDAC substrate** (20 μ M) + 5 μ l BSA (1 mg/ml) + 30 μ l **HDAC11 Assay Buffer**). Add 40 μ l of master mixture to all wells.
- 2) Add 5 μ l of inhibitor solution of each well designated "Test Inhibitor." For the "Positive Control" and "Blank," add 5 μ l of the same solution without inhibitor (inhibitor buffer). Add 5 μ l of diluted **Trichostatin A** (100 μ M) to the wells designated "Inhibitor Control." Keep final DMSO concentration at or below 1%.
- 3) Add 5 μ l of **HDAC11 Assay Buffer** to the wells designated "Blank."
- 4) Initiate reaction by adding 5 μ l 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 μ M)	5 μ l	5 μ l	5 μ l	5 μ l
BSA (1 mg/ml)	5 μ l	5 μ l	5 μ l	5 μ l
HDAC11 Assay Buffer	35 μ l	30 μ l	30 μ l	30 μ l
Diluted Trichostatin A (100 μ M)	-	-	-	5 μ l
Test Inhibitor	-	-	5 μ l	-
Inhibitor buffer (no inhibitor)	5 μ l	5 μ l	-	-
Diluted HDAC11 (10 ng/ μ l)	-	5 μ l	5 μ l	5 μ l
Total	50 μl	50 μl	50 μl	50 μl

Step 2:

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

Step 3:

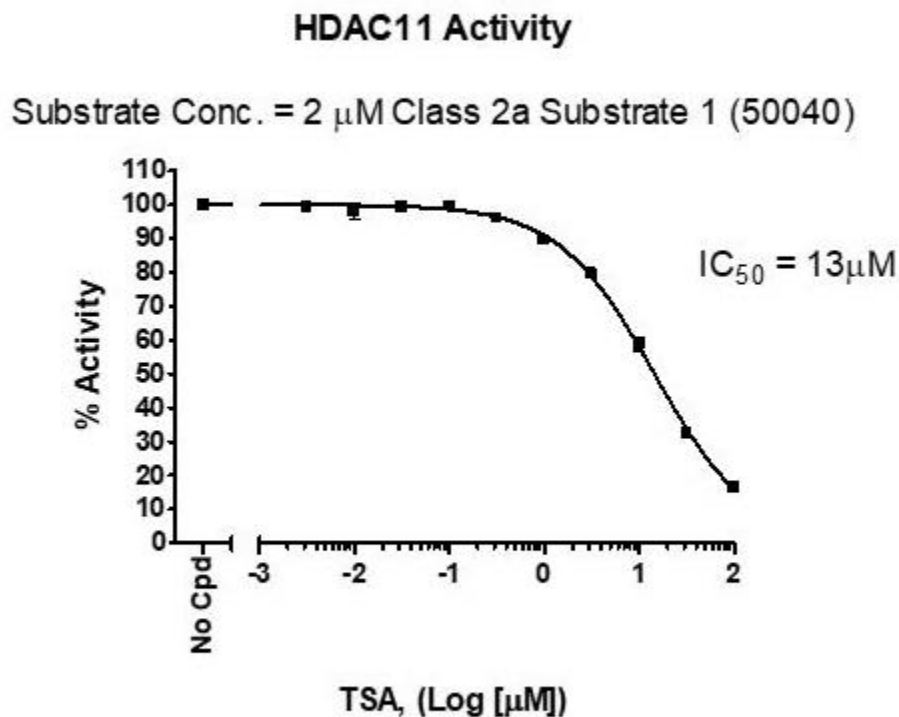
Read sample in a microtiter plate-reading fluorimeter capable of excitation at a wavelength in the range of 350-380 nm and detection of emitted light in the range of 440-460 nm. "Blank" value is subtracted from all other values.

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

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.*

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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
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