



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Kinetic HDAC3 Assay Kit

Catalog #: 53003

DESCRIPTION: The *Kinetic HDAC3 Assay Kit* is a complete assay system designed to measure histone deacetylase 3 (HDAC3) activity kinetically for screening and profiling applications. It comes in a convenient 96-well format, with all the reagents necessary for 100 fluorescent HDAC3 activity measurements. In addition, the kit includes purified HDAC3/NCOR2 enzyme complex and a potent HDAC inhibitor, SAHA, for use as a positive control. The *Kinetic HDAC3 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 one simple step for setting up the HDAC3 reaction on a microtiter plate is needed and then the HDAC3 activity level can be measured kinetically using a fluorescence reader.

COMPONENTS:

Catalog #	Reagent	Amount	Storage	
50003	HDAC3/NCOR2 human recombinant enzyme	2 µg	-80°C	Avoid Freeze/Thaw Cycles!
50037	Fluorogenic HDAC substrate 3 (5 mM)	50 µl	-80°C	
53029	10x HDAC Developer	1 ml	-80°C	
	Developer Dilution Buffer	10 ml	-20°C	
	SAHA (1 mM)	20 µl	-20°C	
50031	HDAC Assay Buffer	20 ml	-20°C	
	BSA (1mg/ml)	1 ml	-20°C	
79685	black, low binding NUNC black microtiter plate	1 plate	Room temp.	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Fluorescent microplate reader

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(S):

1. Santo, L., *et al.*, *Blood*. 2012 Mar 15;**119(11)**:2579-89.
2. Bradner, J.E., *et al.*, *Nat Chem Biol*. 2010 Mar;**6(3)**: 238-243.

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

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

Thaw all of the components except the HDAC enzyme, and place on ice. All of the reactions should be set up on ice.

1) Prepare the master **HDAC Substrate Solution**

To make the master **HDAC Substrate Solution**, mix 2368 μ l **HDAC Assay Buffer**, 800 μ l BSA (1mg/ml), and 32 μ l **HDAC Substrate**. This is enough for 100 reactions. If you have less than 100 reactions, you can reduce the components proportionally. (Make sufficient quantity needed for the assay; store remaining BSA and 5 mM stock solution in aliquots at -80°C.)

Add 20 μ l of the master **HDAC Substrate Solution** to each well.

2) Prepare **Inhibitor Solutions**

Dilute the Test Inhibitor in 10% DMSO at 10-fold higher than the final concentration you want to test (the final DMSO concentration is 1% in all of the reactions).

Dilute the same buffer without the inhibitor (Inhibitor Buffer) in 10% DMSO as a control.

Add 10 μ l of the **Test Inhibitor Solution** to each well designated as "Test Inhibitor", and 10 μ l of **SAHA Solution** to each well designated as "SAHA Inhibitor Control". Add 10 μ l of the Inhibitor Buffer to each well designated "Blank" and "Positive Control".

3) Prepare **Developer Solution**

Dilute the **10X Developer Solution** (1:10 dilution) in **Developer Dilution Buffer**.

Add 50 μ l of the **Developer Solution** to each well.

4) Prepare **HDAC3/NCOR2 Enzyme Solution**

Dilute **HDAC3/NCOR2** in HDAC assay buffer to 0.8 ng/ μ l (16 ng/reaction). Aliquot any remaining enzyme and store undiluted at -80°C.

Add 20 μ l of HDAC assay buffer to each well designed "Blank". Add 20 μ l of the **HDAC3/NCOR2 enzyme solution** to each well designed "Positive Control", "Test Inhibitor" and "SAHA Inhibitor Control". **Always add the enzyme solution last.**

5) Measure the plate in 5 minute intervals for a period of up to one hour, using a microtiter plate-reading fluorimeter capable of excitation in the range of 350-380

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

Please visit our website at: www.bpsbioscience.com

nm and detection of emitted light in the range of 440-460 nm. The “Blank” value is subtracted from all other values.

	“Blank”	Positive Control	Test Inhibitor	SAHA Inhibitor Control
Master HDAC Substrate Solution	20 μ l	20 μ l	20 μ l	20 μ l
SAHA	-	-	-	10 μ l
Test Inhibitor	-	-	10 μ l	-
Inhibitor buffer (no inhibitor)	10 μ l	10 μ l	-	-
1X Developer	50 μ l	50 μ l	50 μ l	50 μ l
HDAC assay buffer	20 μ l	-	-	-
HDAC3/NCOR2 (0.8 ng/ μ l)	-	20 μ l	20 μ l	20 μ l
Total	100 μl	100 μl	100 μl	100 μl

Example of Assay Results:

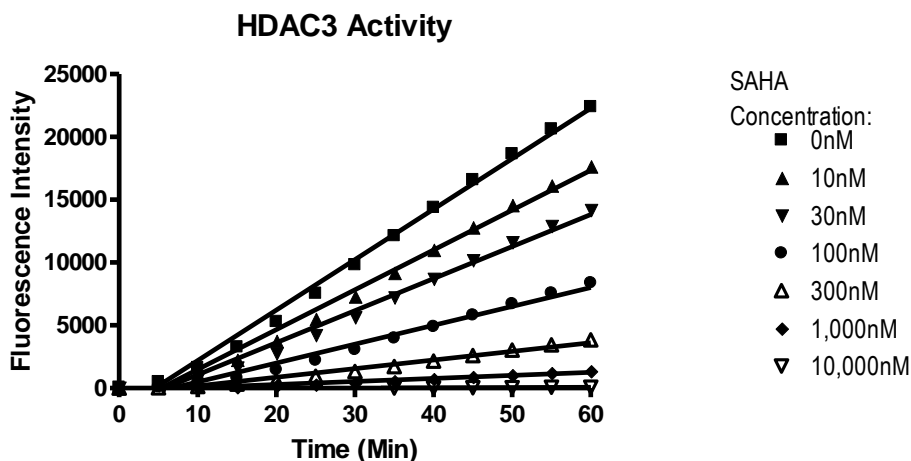


Figure 1. Time course of HDAC3 activity against different concentration of SAH. HDAC3 activity is measured using the *Kinetic HDAC3 Assay Kit*, BPS Bioscience Cat# 53003. Fluorescence was measured using a Tecan fluorescent microplate reader. *Data shown is lot-specific. For lot-specific information, 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

Please visit our website at: www.bpsbioscience.com

HDAC3/NCOR2 Activity

Substrate Conc. = 10 μ M Substrate 3 (50037)

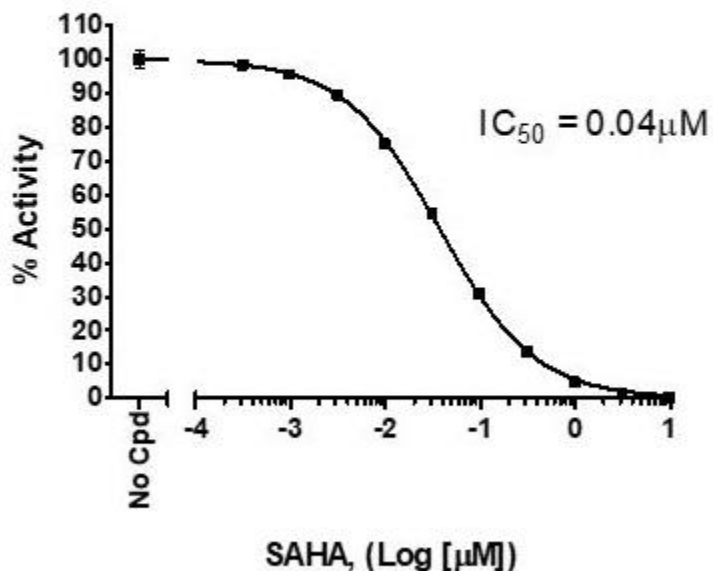


Figure 2. IC₅₀ assay for HDAC3 against SAHA.

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

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
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-Flag)	50052	50 µg
HDAC3/NcoR2	50003	50 µg
HDAC4	50004	10 µg
HDAC5	50005	10 µg
HDAC6 (C-Flag)	50056	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
HDAC1 Assay Kit	50061	96 reactions
HDAC2 Assay Kit	50062	96 reactions
HDAC3 Assay Kit	50073	96 reactions
HDAC4 Assay Kit	50064	96 reactions
HDAC5 Assay Kit	50065	96 reactions
HDAC6 Assay Kit	50076	96 reactions
HDAC7 Assay Kit	50067	96 reactions
HDAC8 Assay Kit	50068	96 reactions
HDAC9 Assay Kit	50069	96 reactions

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

Please visit our website at: www.bpsbioscience.com