



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

Data Sheet

Fluorogenic HDAC/SIRT Substrate

Catalog #: 50032

DESCRIPTION: A fluorogenic, acetylated peptide substrate for HDACs and SIRTs. Based on residues 379-382 of p53 (Arg-His-Lys-Lys(Ac), a site of regulatory acetylation by the p300 and CBP acetyltransferases (lysines 381, 382)¹⁻³, it is a suitable substrate for HDAC2, HDAC3, HDAC6; SIRT1, SIRT2 and SIRT3 tested so far. This substrate should be used in conjunction with HDAC or SIRT Developer, respectively.

PURITY: >90% by HPLC

MOLECULAR WEIGHT: 612.6

APPLICATIONS: Study for enzyme activity, and screening of small molecule inhibitors for drug discovery and HTS applications.

SUPPLIED AS: 5 mM solution in DMSO

QUANTITY: 100 µl

STORAGE: -20°C or -70°C. Protect from exposure to direct light. Avoid freeze/thaw cycles.

REFERENCES:

1. A. Ito *et al.* (2001) *EMBO J.* **20**: 1331.
2. N.A. Barlev *et al.* (2001) *Mol. Cell* **8**: 1243.
3. A. Ito *et al.* (2002) *EMBO J.* **21**: 6236.

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



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

Assay Protocol HDAC:

Materials:

Assay buffer (BPS catalog # 50031)
Assay developer (BPS catalog # 50030)
HDAC Substrate (BPS catalog # 50032)

Step 1: Adding all reaction mixture to a low binding NUNC black plate (VWR catalog number 62408-936)

1 μ l of HDAC2 (0.1 μ g/ μ l)
39 μ l of HDAC assay buffer (BPS catalog # 50031)
5 μ l of 1 mg/ml BSA
5 μ l of 200 μ M substrate (BPS catalog # 50032)

Incubate at 37 °C for 30 min.

Step 2: Stop the reaction

Add 50 μ l of HDAC assay developer (2x) (BPS catalog # 50030) and incubate the plate at room temperature for 15 min.

Step 3:

Read sample in a microtiter-plate reading fluorimeter capable of excitation at a wavelength in the range 350-380 nm and detection of emitted light in the range 440-460 nm.

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



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

Assay Protocol SIRT:

Materials:

SIRT Assay buffer (BPS catalog #50090)
SIRT Assay developer (BPS catalog #50089)
SIRT Substrate (BPS catalog #50080)
SIRT1 (1 µg/µl) (BPS catalog #50012) or other SIRT enzyme

Step 1: Adding all reaction mixture to a low binding NUNC black plate (VWR catalog number 62408-936):

SIRT Assay Buffer, 35 µl
1 mg/ml BSA, 5 µl
200 µM SIRT substrate, 5 µl
SIRT1, 5 µl

Incubate at 37°C for 30 min.

Step 2: Stop the reaction

Add 50 µl of SIRT assay developer (undiluted) and incubate the plate at room temperature for 15 min.

Step 3:

Read samples in a microtiter-plate reading fluorimeter capable of excitation at 350-380 nm and emission at 440-460 nm. "Blank" value (no enzyme negative control) 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
Please visit our website at: www.bpsbioscience.com