

Data Sheet

Fluorogenic HDAC Substrate (Green fluorescence)

Catalog #: 50038 Lot#: 130903

DESCRIPTION: A fluorogenic, acetylated peptide substrate for HDACs. It is a suitable substrate for many HDACs, except class 2a HDACs.

PURITY: >95% by HPLC.

APPLICATIONS: Study enzyme kinetics, and screen small molecule inhibitors for drug discovery and HTS applications.

SUPPLIED AS: DMSO solution

QUANTITY: 50 μl @ 5 mM

STORAGE: -20 ℃ or -70 ℃.

REFERENCES:

- 1. Ito, A., et al., EMBO J. 2001;20:1331.
- 2. Barlev, N.A., et al., Mol Cell. 2001;8:1243.



ASSAY PROTOCOL:

Material: Assay buffer (BPS #50031); Assay developer (BPS #50030); HDAC Substrate (BPS #50038)

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

2 μl of HDAC1 (0.33 μg/μl) 38 μl of HDAC assay buffer (BPS #50031) 5 μl of 1 mg/ml BSA 5 μl of 200 μM substrate (BPS #50038)

Incubate at 37 °C for 30 min.

Step 2: Stop the reaction

Add 50 μ l of HDAC assay developer (2x) (BPS #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 of ~485 nm and detection of emitted light in the range of 528 nm. "Blank" value is subtracted from all other values.