



6044 Cornerstone Court West,
Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email:
info@bpsbioscience.com

Data Sheet

Fluorogenic Sirtuin 5 Substrate

Catalog #: 50126

DESCRIPTION: A fluorogenic, succinylated peptide substrate for Sirtuin 5. This substrate should be used in conjunction with SIRT Developer, #50089. Optimal excitation wavelength 350–380 nm, emission wavelength 440–460 nm.

PURITY: >90% by HPLC.

APPLICATIONS: Suitable for use with BPS's Fluorogenic SIRT5 Assay Kit, #50085, and to study enzyme kinetics and screen small molecule inhibitors of Sirtuin 5 for drug discovery and HTS applications.

SUPPLIED AS: DMSO solution

QUANTITY: 50 µl @ 5 mM

STORAGE: -20°C or -70°C. Avoid freeze/thaw cycles. Stable for at least 6 months from date of receipt when stored as directed. Protect from light.

REFERENCES:

1. Madsen, A.S. and Olsen, C.A., *J. Med. Chem.* 2012, **55**:5582-5590.
2. Du, J., *et al.*, *Science*, 11 November 2011, **334**:806-809.

ASSAY PROTOCOL:

Materials:

SIRT Assay Buffer (BPS # 50090)
SIRT5 Substrate (# 50126)
SIRT5 enzyme (# 50016) or other SIRT5-containing sample
NUNC black 96-well plate, low protein binding (VWR #62408-936).
SIRT Assay Developer (BPS # 50089)

Step 1:

Set up 50 µl of Sirt5 reactions containing SIRT assay buffer, 400 ng SIRT5 enzyme, and 20 µM SIRT5 substrate in the 96-well plate. Incubate at 37°C for 30 min.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** 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 West,
Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email:
info@bpsbioscience.com

Step 2:

Add 50 μ l of SIRT Assay Developer (2x) (BPS #50089) to each well 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. "Blank" value (no enzyme 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 **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com