



Protocol

Assay of HDAC11 (BML-SE560)

with

FLUOR DE LYS[®]-SIRT1 (BML-KI177)

and

FLUOR DE LYS Developer II (BML-KI176)

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

COMPONENTS OF ASSAY:

HDAC Assay Buffer II (BML-KI422)

(50mM Tris/Cl, pH 8.0, 137mM NaCl, 2.7mM KCl, 1mMmgCl₂, 1mg/ml BSA (fatty acid free BSA, e.g. Sigma catalog # A-3803))

HDAC11 (BML-SE560)

Dilute to enough HDAC11 to 0.1µg/µl with HDAC Assay Buffer II (BML-KI422) to provide 5µl per well. Keep on ice until use.

FLUOR DE LYS®-SIRT1 (BML-KI177)

Thaw quickly and keep on ice. Dilute 5mM stock in HDAC Assay Buffer II.

2x Substrate Solution

Prepare a 2x substrate solution by diluting the 5mM “FLUOR DE LYS-**SIRT1**” (BML-KI177) in HDAC Assay Buffer II (BML-KI422). Each assay well will require 25µl (see below). For example, prepare 1ml of 20µM substrate (for final 10µM) by mixing 4µl of 5mM FLUOR DE LYS-**SIRT1** (BML-KI177) and 996µl HDAC Assay Buffer II (BML-KI422). Warm to 37°C before use. (**NOTE:** Determinations of HDAC11’s K_m for “FLUOR DE LYS-**SIRT1**” (BML-KI177) fall in the 5-10µM range. Therefore, a reasonable substrate concentration for inhibitor screening would be 10µM “FLUOR DE LYS-**SIRT1**” (BML-KI177), while 200 or 500µM would be more suitable for a specific activity measurement at a saturating substrate concentration.)

Trichostatin A (BML-GR309; HDAC Inhibitor)

Prepare a 0.2mM stock in dimethylsulfoxide (DMSO). DMSO stock may be stored at -20°C. The 0.2mM stock will be diluted 100-fold in 1x Developer II in order to stop HDAC activity at the start of the signal development process. To prepare a stock for use in testing trichostatin inhibition (i.e. for addition to the deacetylation phase of the reaction), dilute 0.2mM stock to 10µM in HDAC Assay Buffer II (e.g. 5µl plus 95µl) and keep on ice. Addition of 2.5µl of this 10µM stock per well will result in strong inhibition (final [trichostatin A] = 500 nM)

FLUOR DE LYS Developer II (5x Concentrate, Cat. # BML-KI176)

Shortly before use, dilute 5x stock solution to 1x plus 2µM trichostatin A. For example, prepare 1ml by mixing 200µl of the 5x Concentrate, 790µl HDAC Assay Buffer II (BML-KI422) and 10µl 0.2mM trichostatin A in DMSO. Store the 1x Developer II plus trichostatin on ice until use. Do not store excess, but prepare freshly as needed.



Handle
with care

FLUOR DE LYS[®] Deacetylated Standard (BML-KI142)

Dilute the 10mM stock in DMSO to 1 μ M with HDAC Assay Buffer II (BML-KI422).

½ Volume 96-well white NBS micro-plate (BML-KI571)

REACTION CONDITION EXAMPLES :

1. Designate wells for four reactions: 30 min rxn; 30 min rxn plus trichostatin A; 0 min rxn and a Standard well.
2. Add 20 μ l of HDAC Assay Buffer II to the 30 min rxn well and the 0 min rxn well. To the third well (30 min. plus trichostatin) add 2.5 μ l of 10 μ M trichostatin A plus 17.5 μ l of HDAC Assay Buffer II. Allow to equilibrate to assay temperature (37°C). (Leave Standard well empty until step 7).
3. Add 5 μ l of diluted HDAC11 (BML-SE560, 0.1 μ g/ μ l) to the wells for the 0 min, 30 min, and 30 min. plus trichostatin rxns.
4. To start reactions, add 25 μ l of the 2x Substrates (37°C) to both 30 min reaction wells. Allow reactions to run 30 min @ 37°C.
5. Add 50 μ l of 1x Developer II plus trichostatin A to both 30 min. reaction wells.
6. To the 0 min rxn well, add 50 μ l of 1x Developer II plus trichostatin A, immediately followed by 25 μ l of 2x substrate.
7. For the standard, mix 50 μ l of 1 μ M standard with 50 μ l of 1x Developer II plus trichostatin A in the fourth well.
8. Allow 45 min. at 37°C for signal to develop and then read plate in a microplate-reading fluorimeter capable of excitation at a wavelength in the range of 350-380 and detection of emitted light in the range of 440-460 nm.
9. Data analysis: Determine the Δ AFU (Arbitrary Fluorescence Units) for the two 30 min rxns. (AFU of 30 min rxn. (with or without trichostatin) minus AFU of the 0 min rxn). Determine AFU/pmol by dividing the Deacetylated standard reading (AFU) by 50 pmol. Calculate pmol of substrate deacetylated in 30 min (divide Δ AFU by AFU/pmol).

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com