

# FLUOR DE LYS®

## HDAC fluorometric cellular activity assay kit

**Monitor HDAC activity in cell culture using cell-permeable FLUOR DE LYS® substrate**

The HDAC fluorometric cellular activity assay allows the determination of deacetylase activity within an undisturbed cellular environment and provides accurate activity information that is reflective of endogenous regulation. It also allows the study of the effects of upstream regulators on deacetylase activity and the detection of inhibitors or activators that act indirectly to affect deacetylase activity. The FLUOR DE LYS® HDAC substrate is cell permeable. It is deacetylated by cellular HDACs and thus suitable for high-throughput cell-based deacetylase assays. This kit provides the reagents and protocols needed for determining rates of intracellular deacetylase activity with cultured cells. Deacetylation can be quantitated by addition of developer to the media and lysed cells.

HDACs are typically found in multiprotein complexes and are tightly regulated by subcellular localization, phosphorylation, and likely by other mechanisms.

Citations: 19

[View Online »](#)

### Ordering Information

[Order Online »](#)

BML-AK503-0001	96 wells
----------------	----------

### Manuals, SDS & CofA

[View Online »](#)

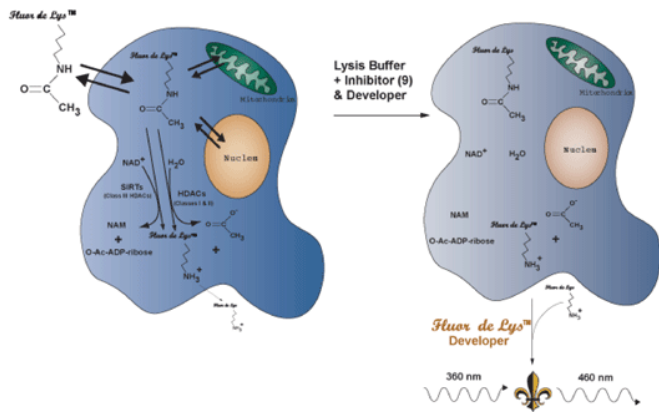


Figure: Method for Assaying Intracellular HDAC Activity with the Fluorogenic, Cell-permeable Substrate, FLUOR DE LYS<sup>®</sup> (Prod. No. BML-KI104). FLUOR DE LYS<sup>®</sup> is added to cell growth medium and enters the cells. Deacetylation by class I/II HDACs or sirtuins yields the deacetylated form of FLUOR DE LYS<sup>®</sup>, some of which may exit the cell. Lysis of the cells with detergent allows contact between the non-cell permeable Developer (Prod. No. BML-KI105) and both intra- and extracellular deacetylated substrate, thus producing a fluorescent signal (symbol). A strong class I/II HDAC inhibitor (e.g. trichostatin A) is added along with the lysis buffer containing detergent to insure that no deacetylation occurs after cell lysis.

# Handling & Storage

**Use/Stability** Store all components except the microplate at -80°C for the highest stability. The HeLa Nuclear Extract, Prod. No. BML-KI345, must be handled with particular care in order to retain maximum enzymatic activity. Defrost it quickly in a RT water bath or by rubbing between fingers, then immediately store on an ice bath. The remaining unused extract should be refrozen quickly, by placing at -80°C. If possible, snap freeze in liquid nitrogen or a dry ice/ethanol bath. To minimize the number of freeze/thaw cycles, aliquot the extract into separate tubes and store at -80°C. The FLUOR DE LYS® Substrate, Prod. No. BML-KI104, when diluted in Assay Buffer, may precipitate after freezing and thawing. It is best, therefore, to dilute only the amount needed to perform the assays of that day.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Cellular Histone deacetylase fluorescent assay kit

**Application** Activity assay, Cell-based assays, Fluorescent detection, HTS

**Nuclear Extracts from Hela Cells** (human cervical cancer cell line) (Prod. No. BML-KI345)

(50 µl; 2 mg protein/ml in 0.1M KCl, 20mM HEPES/NaOH, pH 7.9, 20% (v/v) glycerol, 0.2mM EDTAacid, 0.5M DTT, 0.1mM PMSF; Prepared according to a modification of J.D. Dignam et al (1983) and S.M. Abmayr et al. (1988))

Storage: -80°C, avoid freeze/thaw cycles!

**FLUOR DE LYS<sup>®</sup> Substrate** (Prod. No. BML-KI104)

(50 µl; 50mM in DMSO)

Storage: -80°C

**FLUOR DE LYS<sup>®</sup> Developer Concentrate** (20x) (Prod. No. BML-KI105)

(300 µl; 20x stock solution, dilute in assay buffer before use)

Storage: -80°C

**Trichostatin A** (HDAC Inhibitor) (Prod. No. BML-GR-309-9090)

(100 µl; 0.2mM in DMSO)

Storage: -80°C

**Nicotinamide** (Sirtuin Inhibitor) (Prod. No. BML-KI283)

(500 µl; 50mM nicotinamide in 50mM TRIS/Cl, pH 8.0, 137mM NaCl, 2.7mM KCl, 1mM MgCl<sub>2</sub>)

Storage: -80°C

**FLUOR DE LYS<sup>®</sup> Deacetylated Standard** (Prod. No. BML-KI142)

(30 µl; 10mM in DMSO)

Storage: -80°C

**HDAC Assay Buffer** (Prod. No. BML-KI143)

(50mM TRIS/Cl, pH 8.0, 137mM NaCl, 2.7 mM KCl, 1mM MgCl<sub>2</sub>)

(20 ml)

Storage: -20°C

**Cell Lysis Buffer** (Prod. No. BML-KI346)

(20 ml; 1.0% NP-40 in HDAC Assay Buffer)

Storage: -80°C

**1/2 volume clear microplate** (Prod. No. 80-2404)

Storage: Ambient

**1/2 volume clear microplate, sterile** (Prod. No. 80-2430)

Storage: Ambient

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)