

FLUOR DE LYS®- Green HDAC fluorometric activity assay kit

An improved FLUOR DE LYS® HDAC assay with FLUOR DE LYS®-Green, a new substrate offering higher sensitivity and an excitation and emission (485/530nm) that avoids quenching and fluorescent interference from compounds absorbing in the near UV and blue range. The FLUOR DE LYS®-Green HDAC assay is a complete kit for measuring histone deacetylase (HDAC) activity in cell or nuclear extracts, immunoprecipitates or purified enzymes. It comes in a convenient 96-well format, with all reagents necessary for fluorescent HDAC or sirtuin activity measurements and calibration of the assay. The FLUOR DE LYS®-Green HDAC assay is based on the FLUOR DE LYS®-Green substrate and FLUOR DE LYS® developer combination. The FLUOR DE LYS® system (Fluorogenic Histone deAcetylase Lysyl Substrate/Developer) is a highly sensitive and convenient alternative to radiolabeled, acetylated histones or peptide/HPLC methods for the assay of histone deacetylases. The assay procedure has two steps. First, the FLUOR DE LYS®-Green substrate, which comprises an acetylated lysine side chain, is incubated with a sample containing HDAC activity (HeLa nuclear or other extract, purified enzyme, bead-bound immunocomplex, etc.). Deacetylation of the substrate sensitizes the substrate so that, in the second step, treatment with the FLUOR DE LYS® developer produces a fluorophore.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AK530-0001

96 wells

Manuals, SDS & CofA

[View Online »](#)

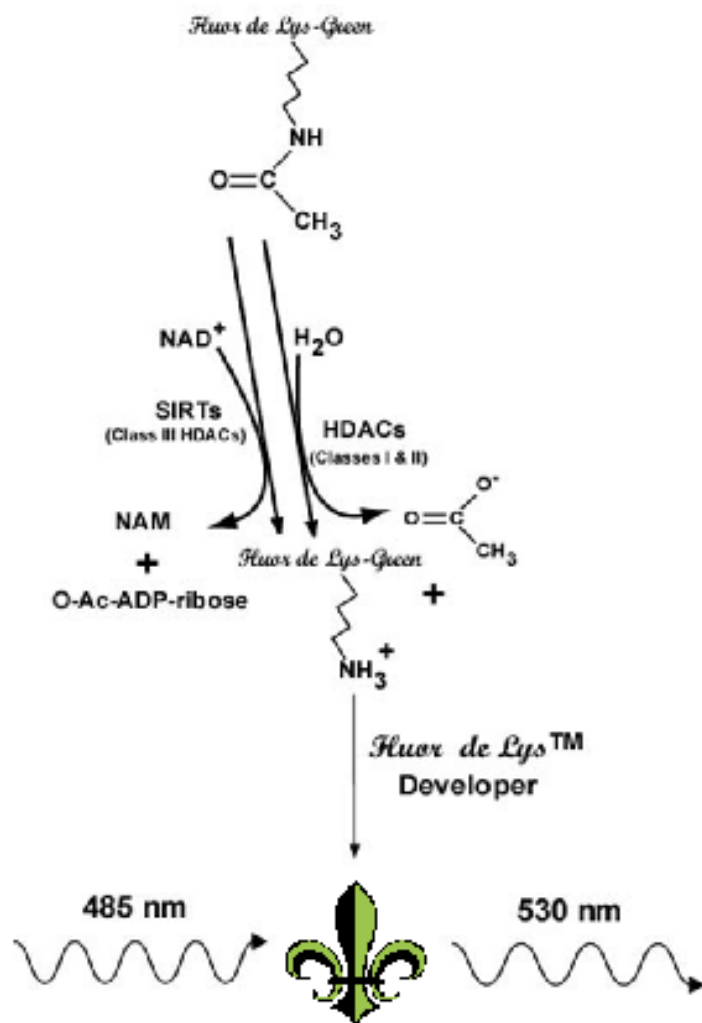


Figure 1: Reaction scheme of the FLUOR DE LYS[®]-Green HDAC assay kit. Deacetylation of the substrate sensitizes it to the developer, which then generates a fluorophore (lily-symbol).

Handling & Storage

Handling	Store all components (except microtiter plates) at -70°C. Unused HeLa Nuclear Extract should be refrozen quickly, if possible snap freeze in liquid nitrogen.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Histone deacetylase green fluorescent assay kit
Application	Activity assay, Cell-based assays, Fluorescent detection, HTS

Contents

Nuclear Extract from HeLa Cells (human cervical cancer cell line) (Prod. No. BML-KI140)

(100µl in 0.1M potassium chloride, 20mM HEPES/NaOH, 20% (v/v) glycerol, 0.2mM EDTA, 0.5mM DTT, 0.5mM PMSF)

Storage: -70°C.

FLUOR DE LYS[®]-Green Substrate (Prod. No. BML-KI572)

(50µl 50mM in DMSO)

Storage: -70°C.

FLUOR DE LYS[®] Developer Concentrate (20X) (Prod. No. BML-KI105)

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

Storage: -70°C.

Trichostatin A (Prod. No. BML-GR309)

(100µl 0.2mM in DMSO)

Storage: -70°C.

FLUOR DE LYS[®]-Green Standard (Prod. No. BML-KI605)

(30µl 1mM in DMSO)

Storage: -70°C.

Note: BML-KI605 is an improved replacement of BML-KI573.

NAD⁺ (Sirtuin Substrate) (Prod. No. BML-KI282)

(500µl; 50mM in 50mM TRIS, pH 8.0, 137mM sodium chloride, 2.7mM potassium chloride, 1mM magnesium chloride)

Storage: -70°C.

Nicotinamide (Prod. No. BML-KI283)

(500µl; 50mM in 50mM TRIS, pH 8.0, 137mM sodium chloride, 2.7mM potassium chloride, 1mM magnesium chloride)

Storage: -70°C.

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

(20ml; 50mM TRIS, pH 8.0, 137mM sodium chloride, 2.7mM potassium chloride, 1mM magnesium chloride)

Storage: -70°C.

96 Well Plates (Prod. No. BML-KI101)

Storage: Room temperature

White NBS Microplate (Prod. No. BML-KI571)

Storage: Room temperature

Technical Info / Product Notes

Biocompare Article: [Keep an Eye on Apoptosis with Caspase Assays](#)



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-
fr@enzolifesciences.com](mailto:info-fr@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)