

# FLUOR DE LYS<sup>®</sup>

## developer I

## concentrate

### An essential reagent for signal generation in HDAC or sirtuin assays done with the FLUOR DE LYS<sup>®</sup> substrate.

FLUOR DE LYS<sup>®</sup> Developer is an essential reagent for signal generation in HDAC or sirtuin assays done with the FLUOR DE LYS<sup>®</sup> substrate. An appropriate FLUOR DE LYS<sup>®</sup> family substrate, comprising an acetylated lysine side chain, is first incubated with a sample containing HDAC activity (nuclear or cellular extract, purified enzyme, bead-bound immunocomplex, etc.). Deacetylation of the substrate sensitizes it so that, in a second step, treatment with the FLUOR DE LYS<sup>®</sup> Developer produces a fluorophore. Typically, an 'inhibitor stop' will be included in the 1x Developer dilution (e.g. Trichostatin A) for class I and II HDACs, nicotinamide for sirtuins).

Citations: 8

[View Online »](#)

### Ordering Information

[Order Online »](#)

BML-KI105-0300	300µl
----------------	-------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Concentration** 20x concentrate

**Quantity** After dilution (1x), sufficient for 100 assays (50 µl reaction volume).

**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)