

# FLUOR DE LYS<sup>®</sup>

## developer II

## concentrate

FLUOR DE LYS<sup>®</sup> Developer II is an essential reagent for signal generation in assays done with the FLUOR DE LYS<sup>®</sup>-HDAC8, FLUOR DE LYS<sup>®</sup>-SIRT1 and FLUOR DE LYS<sup>®</sup>-SIRT2 substrates. Deacetylation of the FLUOR DE LYS<sup>®</sup> substrate sensitizes it so that, in a second step, treatment with the FLUOR DE LYS<sup>®</sup> Developer II produces a fluorophore. Typically, an 'inhibitor stop' will be included in the 1x Developer dilution (e.g. nicotinamide for sirtuins).

Citations: 23

[View Online »](#)

### Ordering Information

[Order Online »](#)

BML-KI176-1250	1.25ml
----------------	--------

### 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

### Quantity

NOTE: This product is supplied packaged in 5 tubes (250 µl per tube) for a total volume of 1.25 ml. This packaging is used for optimal stability of the developer.

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)