

FLUOR DE LYS[®]

developer I

concentrate

An essential reagent for signal generation in HDAC or sirtuin assays done with the FLUOR DE LYS[®] substrate.

FLUOR DE LYS[®] Developer is an essential reagent for signal generation in HDAC or sirtuin assays done with the FLUOR DE LYS[®] substrate. An appropriate FLUOR DE LYS[®] 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[®] 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).



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)