

FLUOR DE LYS[®]

deacetylase substrate

Highly sensitive and convenient alternative to radiolabeled, acetylated histones or peptide/HPLC methods for the assay of histone deacetylases.

FLUOR DE LYS[®] Substrate is a highly sensitive and convenient alternative to radiolabeled, acetylated histones or peptide/HPLC methods for the assay of histone deacetylases. The FLUOR DE LYS[®] Substrate, which comprises an acetylated lysine side chain, is incubated with a sample containing HDAC activity (nuclear or cellular extract, purified enzyme, bead-bound immunocomplex, etc.). Deacetylation of the substrate sensitizes the substrate so that, in a second step, treatment with the FLUOR DE LYS[®] Developer produces a fluorophore. Histone deacetylases from all known classes have been successfully assayed with the FLUOR DE LYS[®] Substrate. These include human HDACs 1, 2, 3 and 8 (class I), human HDACs 4 and 9 (class II) and the sirtuins yeast Sir2, human SIRT1 and human SIRT2 (class III; reaction must include the sirtuin substrate NAD⁺). The substrate is cell-permeable and serves as the basis for the HDAC Cellular Activity Assay Kit (Prod. No. BML-AK503). **NOTE:** Must be used in conjunction with FLUOR DE LYS[®] Developer (Prod. No. BML-KI105). Sufficient for 100-1000 assays of HDACs or sirtuins (50 µl reaction volume, 50-500 µM (substrate)).

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-KI104-0050	50µl
----------------	------

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

Formulation Supplied as a 50 mM solution (50 µl) in dimethylsulfoxide (DMSO).

MW 387.4

Purity ≥95% (HPLC)

Quantity 2.5 µmol (50 µl)



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)