

FLUOR DE LYS[®] -

Green substrate

The FLUOR DE LYS[®]-Green substrate offers 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 (Prod. No. BML-KI105) produces a fluorophore. The FLUOR DE LYS[®]-Green substrate has a higher wavelength excitation and emission than the original FLUOR DE LYS[®] substrate (Prod. No. BML-KI104). This avoids interference by quenching or fluorescence from compounds absorbing and/or emitting in the near UV and blue range.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-KI572-0050	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

Couple Target HDAC

Couple Type Substrate

Formulation 50mM solution in DMSO.

Purity ≥95% (HPLC)

Quantity 2.5 µmol

Technical Info / Product Notes **NOTE:** Must be used in conjunction with FLUOR DE LYS[®] Developer (Prod. No. BML-KI105; see attached assay conditions).



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)