

FLUOR DE LYS[®]- HDAC8 deacetylase substrate

A fluorogenic, diacetylated peptide substrate for HDAC8 (histone deacetylase-8).

FLUOR DE LYS[®]-HDAC8 is a fluorogenic, diacetylated peptide substrate for HDAC8 (histone deacetylase-8). Based on residues 379-382 of p53 (Arg-His-Lys(Ac)-Lys(Ac)), a site of regulatory acetylation by the p300 and CBP acetyltransferases (lysines 381, 382), it was the best for HDAC8 from among a panel of substrates patterned on p53, histone H3 and histone H4 acetylation sites. FLUOR DE LYS[®]-HDAC8 is deacetylated by HDAC8 at a rate of more than 10-fold that of the acetylated lysine substrate, FLUOR DE LYS[®] (BML-KI104; substrates both at 100 μ M). Although named because of HDAC8's preference for it, it is also an excellent substrate for SIRT1 (BML-SE239; 5x the rate of BML-KI104 at 25 μ M, 0.5 mM NAD⁺), SIRT2 and HeLa Nuclear Extract (BML-KI140; 3x KI-104 rate, 25 μ M). Must be used in conjunction with FLUOR DE LYS[®] Developer II (BML-KI176). Sufficient for 100-200 assays of human recombinant HDAC8 (BML-SE145; 1 U/well, 50-100 μ M substrate).

Citations: 19

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-KI178-0005	0.5 μ mol
----------------	---------------

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 5 mM solution (100 µl) in HDAC Assay Buffer.

Purity ≥95%



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)