

FLUOR DE LYS[®]-SIRT2 deacetylase substrate

FLUOR DE LYS[®]-SIRT2 is a fluorogenic peptide substrate for SIRT2 (human Sirtuin 2). Based on residues 317-320 of p53 (Gln-Pro-Lys-Lys(Ac)), a site of regulatory acetylation by the PCAF acetyltransferase (lysine 320), it was the best for SIRT2 from among a panel of substrates patterned on p53, Histone H3, and Histone H4 acetylation sites. FLUOR DE LYS[®]-SIRT2 deacetylase substrate is deacetylated by SIRT2 at a rate of more than 50-fold that of the acetylated lysine substrate (Prod. No. BML-KI104; acetylated substrates both at 25 μ M, 500 μ M NAD⁺). Must be used in conjunction with FLUOR DE LYS[®] Developer II (Prod. No. BML-KI176). Sufficient for 100 assays of human recombinant SIRT2 (Prod. No. BML-SE251; 2U/well, 100 μ M substrate).

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-KI179-0005	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 5mM solution (100 µl) in HDAC Assay Buffer.

Purity ≥95% (HPLC)



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)