

# FLUOR DE LYS®-SIRT1 deacetylase substrate

FLUOR DE LYS®-SIRT1 is a fluorogenic, acetylated peptide substrate for SIRT1 (human Sirtuin 1). Based on residues 379-382 of p53 (Arg-His-Lys-Lys(Ac)), a site of regulatory acetylation by the p300 and CBP acetyltransferases (lysines 381, 382), it was the best substrate for SIRT1 from among a panel of substrates patterned on p53, histone H3, and histone H4 acetylation sites. FLUOR DE LYS®-SIRT1 is deacetylated by SIRT1 (BML-SE239) at a rate of more than 8-fold that of the acetylated lysine substrate, FLUOR DE LYS® (Prod. No. BML-KI104; acetylated substrates both at 25  $\mu$ M, 500  $\mu$ M NAD<sup>+</sup>). The  $K_m$  of FLUOR DE LYS®-SIRT1 for human recombinant Sirtuin 1 is 108  $\mu$ M (determined at 37°C, 500  $\mu$ M NAD<sup>+</sup>). Must be used in conjunction with FLUOR DE LYS® Developer II (Prod. No. BML-KI176). Fluorescent signal indicates deacetylation of Lys382. Sufficient for 100-200 assays of human recombinant SIRT1 (1U/well, 50-100  $\mu$ M substrate).

Citations: 14

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-KI177-0005	0.5 $\mu$ 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 in HDAC Assay Buffer.

**Purity** ≥95% (HPLC)

**Quantity** 0.5 µmol (100 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)