

# FLUOR DE LYS® -H4-AcK16 deacetylase substrate

The deacetylated state of lysine-16 of histone H4 (H4-K16) is thought to be a critical factor in determining regions of chromatin silencing. Conversely, acetylation of lysine-16 can be associated with increased transcription, such as that due to dosage compensation from the X-chromosome of male *Drosophila*. The acetylated H4 lysine-16 is preferentially deacetylated by yeast Sir2 *in vitro* and *in vivo*. FLUOR DE LYS®-H4-AcK16 is a fluorogenic, acetylated peptide substrate suitable for various lysine deacetylases, including the HDAC activities in HeLa nuclear extract, HDAC8, SIRT1 (human Sirtuin 1) and SIRT2 (human Sirtuin 2). It is based on residues from histone H4 (12-16, Lys-Gly-Gly-Ala-Lys(Ac)) and was the preferred substrate for both SIRT1 and SIRT2 within a group patterned on acetylation sites in histone H3 and histone H4 N-terminal tails. FLUOR DE LYS®-H4-AcK16 is deacetylated by SIRT1 (BML-SE239) at a rate of more than 6-fold and by SIRT2 (Prod. No. BML-SE251) at a rate of more than 40-fold those achieved with the acetylated lysine substrate, FLUOR DE LYS® (Prod. No. BML-KI104; acetylated substrates both at 25 µM, 500 µM NAD ). Must be used in conjunction with FLUOR DE LYS® Developer (Prod. No. BML-KI105). Sufficient for 100-200 assays (50-100 µM substrate).

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-KI174-0005	0.5µmol
----------------	---------

## 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.

**MW** 298.4

**Purity** ≥95%

**Last modified: May 29, 2024**



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)