

SCREEN-WELL®

Epigenetics library

The SCREEN-WELL® Epigenetics library contains 43 compounds with defined activity against enzymes which carry out epigenetic modification of lysine. It also includes DNA methylation inhibitors. Compounds are dissolved in DMSO at 10mM and aliquoted into deep-well plates at 100 or 500µl per well. The library is a useful tool for chemical genomics, assay development and other pharmacological applications. The library contains inhibitors of these important enzymes: HDACs, SIRT6, Lysine demethylases, HATs, Histone methyl transferases, DNA methyltransferases as well as SIRT6 activators. A variety of structurally and mechanistically different compound classes are included.

Citations: 25

[View Online »](#)

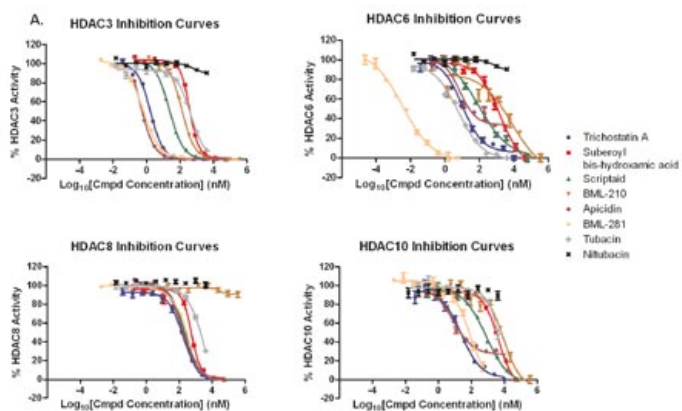
Ordering Information

[Order Online »](#)

BML-2836-0100	100 µl/well - 1Library
BML-2836-0500	500 µl/well - 1Library

Manuals, SDS & CofA

[View Online »](#)



Download our Application Note: A Flexible and Robust Solution for Automated 384-well HDAC Profiling.”””

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year from the date of receipt when stored at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	HTS
Application Notes	Useful tool for chemical genomics, assay development, or other pharmacological applications in the study of epigenetic modifications of lysine and DNA methylation.
Contents	43 compounds supplied in DMSO at a concentration of 10mM.



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)