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, SIRTs, Lysine demethylases, HATs, Histone methyl transferases, DNA methyltransferases as well as SIRT activators. A variety of structurally and mechanistically different compound classes are included.

Citations: 25

View Online »

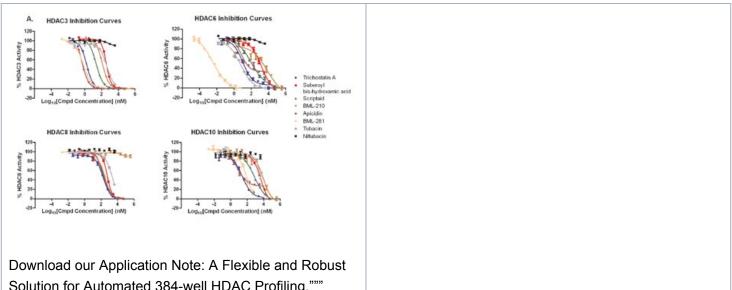
Ordering Information

Order Online »

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

Manuals, SDS & CofA

View Online »



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.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com