

# EPIGENETIC SERVICES

Epigenetics is the study of transmissible modifications in gene expression that occur independently of mutations to the DNA sequence. The deregulation of epigenetic control is a common feature of a number of syndromes, disorders and cancers. This field of research is now considered to be a novel avenue for chemical safety assessment and drug discovery.

## ERASERS

ASSAY	TECHNOLOGY / REF.
DEMETHYLASES	TR-FRET
FBXL10	3695
FBXL11	3596
JARID1A	3697
JARID1B	3701
JARID1C	3703
JARID1D	3696
JMJD1A	3597
JMJD2A	3595
JMJD2B	3598
JMJD2C	3599
JMJD2D	3600
JMJD2E	3601
JMJD3	3602
LSD1	3712

ASSAY	TECHNOLOGY / REF.
PHF8	3702
UTX	3705
<b>HDACS &amp; SIRTUINS</b>	<b>Fluorimetry</b>
HDAC1	2491
HDAC2	2492
HDAC3	2083
HDAC4	2493
HDAC5	2494
HDAC6	2495
HDAC7	2610
HDAC8	2247
HDAC9	2611
HDAC10	2662
HDAC11	2663

ASSAY	TECHNOLOGY / REF.
Sirtuin 1	2991/2581
Sirtuin 2	2582
Sirtuin 3	2583
Sirtuin 6	3808
Sirtuin 7	3809
<b>UBIQUITIN MODIFYING ENZYMES</b>	<b>LanthaScreen™</b>
BAP1	3890
USP 3	3883
USP 7	3887
USP 10	3884
USP 14	3888
USP 16	3889
USP 21	3873

## WRITERS

ASSAY	TECHNOLOGY / REF.
<b>DNA METHYLTRANSFERASES</b>	<b>Radioactivity</b>
DNMT1	3416
DNMT3a	3706
DNMT3b	3708
DNMT3b/3L	3775
<b>KINASES</b>	<b>Radioactivity</b>
Aurora B*	3812
DAPK3/ZIP*	3813
Haspin*	3814
IKKa*	3789
MSK1*	3792
MSK2*	3793
Pim1*	3794
PKBa/Akt1*	3795
PKBb/Akt2*	3796
PKBg/Akt3*	3797
RSK2*	3798
<b>HISTONE ACETYLTRANSFERASES</b>	<b>Radioactivity</b>
CREBBP	3783
EP300	3710
HAT1	3892
KAT2A/GCN5L2	3634

ASSAY	TECHNOLOGY / REF.
KAT5/TIP60	3776
MYST2	3768
MYST3	3766
MYST4	3767
PCAF	3737
<b>HISTONE METHYLTRANSFERASES</b>	<b>Radioactivity</b>
ASH1L	3903
DOT1L	3565
EHMT1	3566
EZH1/EED/SUZ12	3692
EZH2/EED/SUZ12 (PRC2 complex)	3578
G9a	3560
MLL complex	3567
MLL2 complex	3898
MLL3 complex	3899
MLL4 complex	3900
PRDM9	3715
PRMT1	3707
PRMT3	3577
PRMT4	3711
PRMT5 complex	3568
PRMT6	3698
PRMT7	3894

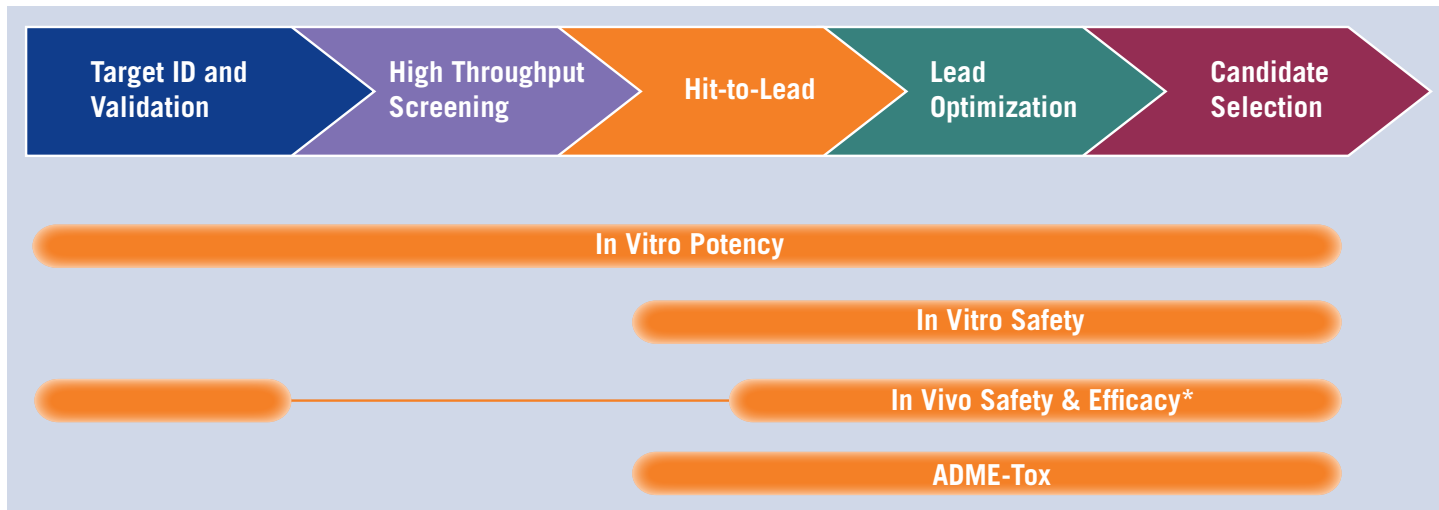
ASSAY	TECHNOLOGY / REF.
PRMT8	3765
SET1B complex	3910
SETD2	3716
SETD7	3569
SETD8	3570
SETDB1	3620
SMYD2	3717
SUV39H1	3704
SUV39H2	3571
SUV420H2	3579
NSD1	3694
WHSC1/NSD2	3749
WHSC1L1/NSD3	3764
<b>SMMTs</b>	<b>Mass Spectrometry</b>
COMT	3946
GAMT	3947
GNMT	3949
HNMT	3943
NNMT	3951
PNMT	3945
TPMT	3942

\* Assays developed using relevant epigenetic substrate.

All assays are human except DNMT3b, ref. 3708 (mouse).

## Comprehensive Drug Discovery Products and Services

In addition to Epigenetic Services, Eurofins Pharma Discovery Services offers a comprehensive package of other drug discovery solutions. This extensive range of products and services aids in the decision making process, starting with Target ID and Validation through to Candidate Selection. Key products and services include:



### In Vitro Potency

Target-based in vitro binding and functional assays for preliminary assessments of efficacy.

### In Vivo Safety & Efficacy

240+ disease-relevant in vivo models to evaluate drug efficacy, along with models for early PK and toxicity assessment through our Partner Lab Pharmacology Discovery Services\*.

\*In vivo Services are conducted by Pharmacology Discovery Services, a Partner Lab of Eurofins Pharma Discovery Services

### In Vitro Safety

Early assessment across a range of targets for safety and toxicity, including: CardiacProfiler™, SafetyScreen44™ and SafetyScreen87™.

### ADME-Tox

140+ ADME-Tox assays to help clients characterize and better understand the DMPK properties of a drug candidate.

For more information visit: [eurofins.com/epds](http://eurofins.com/epds)

Eurofins Panlabs, Inc  
15 Research Park Drive  
St Charles, MO 63304

#### USA

T | +1 (844) 522 7787

F | +1 (636) 362 7131

E | [discoveryservices@eurofins.com](mailto:discoveryservices@eurofins.com)

Eurofins Pharma Discovery  
Services UK, Ltd.  
Gemini Crescent, Dundee Technology Park  
Dundee, DD2 1SW

#### UNITED KINGDOM

T | +44 1382 561600

F | +44 1382 561601

E | [discoveryservices@eurofins.com](mailto:discoveryservices@eurofins.com)

Eurofins Cerep SA  
Le bois l'Évêque  
86600 Celle l'Evescault

#### FRANCE

T | +33 (0)5 49 89 30 00

F | +33 (0)5 49 43 21 70

E | [cerepanlabs@eurofins.com](mailto:cerepanlabs@eurofins.com)

Eurofins Panlabs Taiwan, Ltd.  
No. 158 Li Teh Road, Peitou  
Taipei, 11259

#### TAIWAN

T | +886 2 2898 7888

F | +886 2 2894 8267

E | [discoveryservices@eurofins.com](mailto:discoveryservices@eurofins.com)