

EPIGENETIC PROFILING

The only CRO to bring you validated reader, writer and eraser assays in epigenetics

	Protein name	Reference	Technology
READER	Bromodomain	BAZ2B	AlphaScreen®
	Bromodomain	BRD2(1)	
	Bromodomain	BRD4(1)	
	Bromodomain	CREBBP	
	Bromodomain	FALZ	
ERASER	Demethylase	JMJD1A	LANCE®
	Demethylase	JMJD2A	
	Demethylase	JMJD2C	
	Demethylase	JMJD2E	
	Demethylase	JMJD3	
	HDAC & Sirtuins	HDAC2	Fluorimetry
	HDAC & Sirtuins	HDAC3	
	HDAC & Sirtuins	HDAC4	
	HDAC & Sirtuins	HDAC5	
	HDAC & Sirtuins	HDAC6	
	HDAC & Sirtuins	HDAC7	
	HDAC & Sirtuins	HDAC8	
	HDAC & Sirtuins	HDAC9	
	HDAC & Sirtuins	HDAC10	
	HDAC & Sirtuins	HDAC11	
HDAC & Sirtuins	Sirtuin 1	2991/2581	
HDAC & Sirtuins	Sirtuin 2	2582	
HDAC & Sirtuins	Sirtuin 3	2583	
WRITER	Methyltransferase	DOT1L	Radioactivity
	Methyltransferase	EHMT1	
	Methyltransferase	G9α	
	Methyltransferase	MLL complex	
	Methyltransferase	PRC2 complex	
	Methyltransferase	PRMT3	
	Methyltransferase	PRMT5 complex	
	Methyltransferase	SETD7	
	Methyltransferase	SETD8	
	Methyltransferase	SETDB1 (available soon)	
	Methyltransferase	SUV39H2	

Anticipating
clinical effects
from
in vitro data