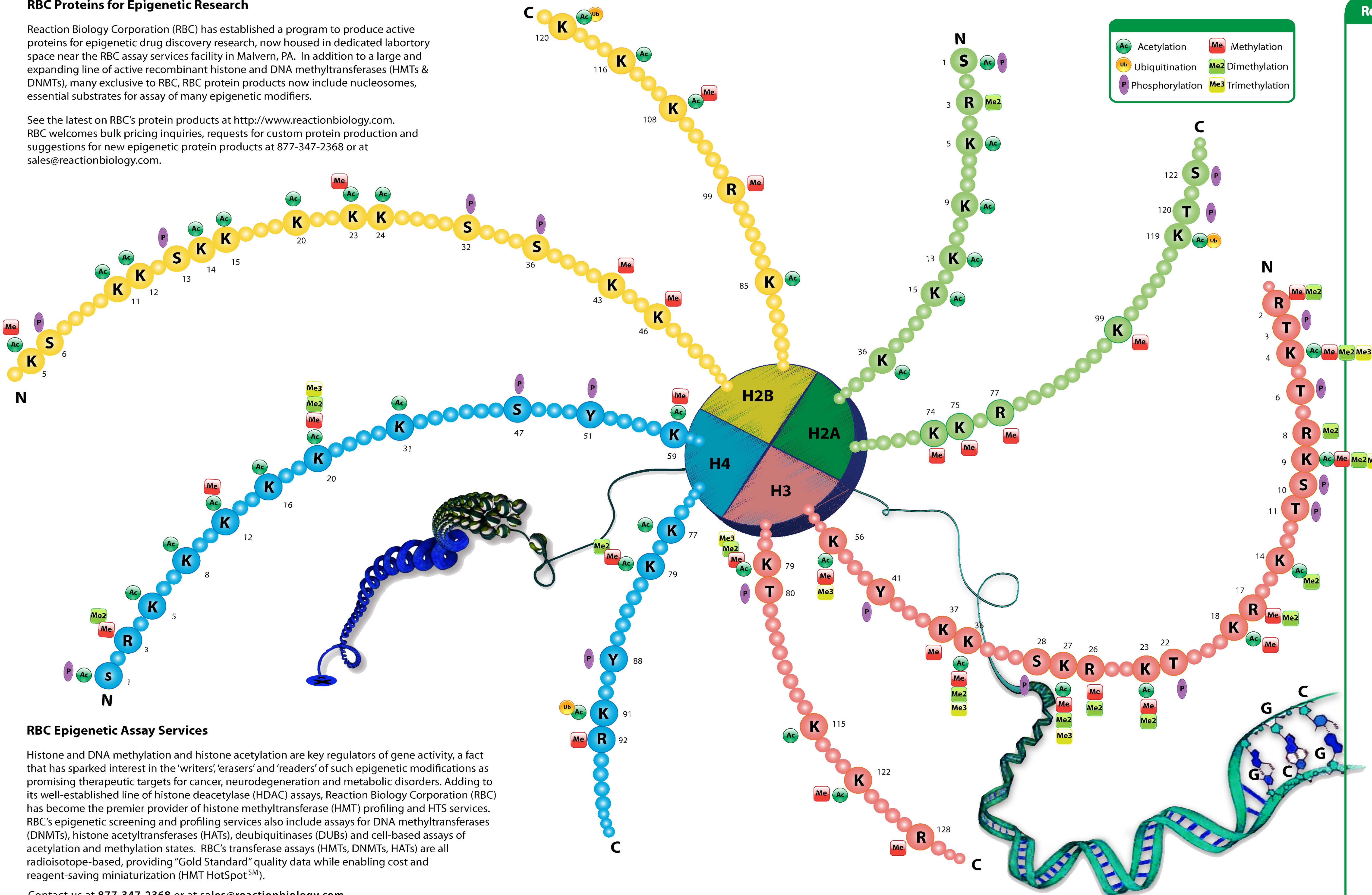


Histone Modification

RBC Proteins for Epigenetic Research

Reaction Biology Corporation (RBC) has established a program to produce active proteins for epigenetic drug discovery research, now housed in dedicated laboratory space near the RBC assay services facility in Malvern, PA. In addition to a large and expanding line of active recombinant histone and DNA methyltransferases (HMTs & DNMTs), many exclusive to RBC, RBC protein products now include nucleosomes, essential substrates for assay of many epigenetic modifiers.

See the latest on RBC's protein products at <http://www.reactionbiology.com>. RBC welcomes bulk pricing inquiries, requests for custom protein production and suggestions for new epigenetic protein products at 877-347-2368 or at sales@reactionbiology.com.



RBC Epigenetic Assay Services

Histone and DNA methylation and histone acetylation are key regulators of gene activity, a fact that has sparked interest in the 'writers,' 'erasers' and 'readers' of such epigenetic modifications as promising therapeutic targets for cancer, neurodegeneration and metabolic disorders. Adding to its well-established line of histone deacetylase (HDAC) assays, Reaction Biology Corporation (RBC) has become the premier provider of histone methyltransferase (HMT) profiling and HTS services. RBC's epigenetic screening and profiling services also include assays for DNA methyltransferases (DNMTs), histone acetyltransferases (HATs), deubiquitinases (DUBs) and cell-based assays of acetylation and methylation states. RBC's transferase assays (HMTs, DNMTs, HATs) are all radioisotope-based, providing "Gold Standard" quality data while enabling cost and reagent-saving miniaturization (HMT HotSpotSM).

Contact us at 877-347-2368 or at sales@reactionbiology.com.

Reaction Biology Epigenetic Assay Services

RBC Assay Services
RBC Proteins Available
Both Assay Services & Proteins Available

DNA Methyltransferases (DNMTs)
DNMT1
DNMT3a
DNMT3b
DNMT3b/3L Complex

Histone (Protein) Methyltransferases (HMTs)

Ash1L	PRDM8
Dot1L	PRDM9
EZH1 Complex	PRMT1
EZH2 Complex	PRMT3
EZH2-Y641F Complex	PRMT4
G9a	PRMT5/MEP50 Complex
GLP	PRMT6
G9a-GLP Complex	PRMT8
MLL1 Complex	SET1A Complex
MLL2 Complex	SET1B Complex
MLL3 Complex	SET7/9
MLL4 Complex	SET8
MLL5 Complex	SETD2 (GST)
NSD1	SETD2 (His)
NSD2	SETMAR
NSD2 (His)	SMYD2
NSD2-E109K	SUV39H1
NSD2-T1150A	SUV39H2
NSD3	SUV420H1-tv1
	SUV420H2-tv2
	WRAD2 Complex

Substrates
GST-GAR
Histone H2A
Histone H2B
Histone H3.3
Nucleosomes (HeLa Mono/Di)
Nucleosomes (HeLa Oligo)
Nucleosomes (HeLa Oligo, Biotinylated)

Histone Deacetylases (HDACs)

HDAC1	HDAC9
HDAC2	HDAC10
HDAC3	HDAC11
HDAC4	Nuclear Extract
HDAC5	SIRT1
HDAC6	SIRT2
HDAC7	SIRT3
HDAC8	SIRT5

Histone Acetyltransferases (HATs)

P300
CBP
hGCN5
KAT5
MYST2
MYST4
pCAF

Reader Domains

ASH1L-[BRD]	BRDT-2	SMARCA2a
ATAD2	BRDT-Tandem	SMARCA2b
ATAD2B	BRPF1a	SMARCA4
BAZ1A	BRPF1b	SP100
BAZ1B	BRPF3	SP110c
BAZ2A	BRWD1-2	SP140c
BAZ2B	CECR2	SP140L
BPTF-[BRD]	CREBBP	TAF1-1
BRD1	EP300	TAF1-2
BRD2-1	HP1β	TAF1L-1
BRD2-2	HP1β-[CHR]	TAF1L-2
BRD2-Tandem	KAT2A	TAF1L-Tandem
BRD3-1	KAT2B	TRIM24
BRD3-2	L3MBTL1	TRIM28
BRD3-Tandem	PB1-1	TRIM33a (His)
BRD4 Full Length	PB1-2	TRIM33b
BRD4-1	PB1-3	TRIM66
BRD4-2	PB1-4	UHRF1-[PHD]
BRD4-Tandem	PB1-5	UHRF1-[SRA]
BRD7	PB1-6	UHRF1-[TDR]
BRD9	PHIP-2	UHRF1-[TDR-PHD]
BRDT-1	PHIP-Tandem	UHRF1-FL

Demethylases

LSD1

Deubiquitinases (DUBs)

A20	UCHL3
Ataxin3	USP10
BAP1	USP14
NEDP1	USP2
SEN1	USP5
UCHL1	USP8

Cellular Assays

H3K27me3
H3K4me2
H3K79me2
H3K36me2
H3K9me2
H3K27me2
Histone Acetylation
Tubulin Acetylation

Other
OGT (O-GlcNAc Transferase)