



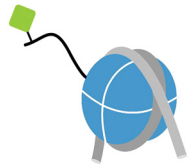
EpiCypher[®]

Bringing Epigenetics to Life

Functionalized
Recombinant Nucleosomes
For Drug Discovery and
Chromatin Research

Functionalized Nucleosome Substrates for Drug Discovery and Chromatin Research

Nucleosomes are the physiological target of readers, writers and erasers that interact with or modify chromatin. The incorporation of nucleosome substrates into drug discovery assays is a dramatic improvement over peptides, providing access to historically challenging targets.



Features

dNuc[™] Designer Nucleosomes

- Contain physiological histone PTMs
- 601 Nucleosome positioning sequence (biotinylated or non-biotinylated)

Benefits

- Suitable for enzyme assays and high-throughput screening
- Suitable for protein-protein interaction studies involving the modification of interest



rNuc[™] Recombinant Nucleosomes

- Fully recombinant human histones
- 601 Nucleosome positioning sequence (biotinylated or non-biotinylated)

- Devoid of post-translational modifications
- Suitable for enzyme assays, inhibitor testing and high-throughput screening



oncoNuc[™] Oncogenic Nucleosomes

- Contains specific mutations associated with cancer
- 601 Nucleosome positioning sequence (biotinylated or non-biotinylated)

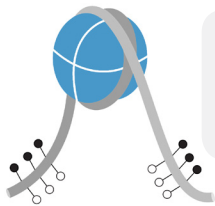
- Study effects of mutations on enzyme activity
- Suitable for high-throughput screening and inhibitor testing



vNuc[™] Histone Variant Nucleosomes

- Includes one of several histone variants
- 601 Nucleosome positioning sequence (biotinylated or non-biotinylated)

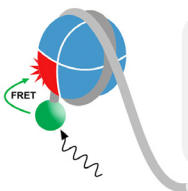
- Study effects of mutations on enzyme activity
- Suitable for high-throughput screening and inhibitor testing



methyl DNA Nuc

- Hemi-methylated 601 Nucleosome positioning sequence
- Methylated DNA can be assembled on modified octamers upon request

- Study effects of DNA methylation on enzyme activity
- Suitable for enzyme assays, inhibitor testing and high-throughput screening



EpiDyne[®] Chromatin Remodeling Assay Substrates

- Nucleosome positioning sequence with an added nucleosome acceptor sequence
- Functionalized DNA or histones to enable HTS assay development

- Suitable for high-throughput screening and inhibitor testing
- Can be used for structural studies



SNAP-ChIP[®] Panels

- DNA-barcoded 601 Nucleosome positioning sequence
- Contain physiological histone PTMs

- Quantitative spike-in controls for ChIP
- Antibody specificity testing
- Monitor experimental variability

dNucs: Designer Recombinant Nucleosomes With PTMs (Biotinylated)*



dNucs

Histone Lysine Methylation

H3K4me1	16-0321	50 µg
H3K4me2	16-0334	50 µg
H3K4me3	16-0316	50 µg
H3K9me1	16-0325	50 µg
H3K9me2	16-0324	50 µg
H3K9me3	16-0315	50 µg
H3K27me1	16-0338	50 µg
H3K27me2	16-0339	50 µg
H3K27me3	16-0317	50 µg
H3K36me1	16-0322	50 µg
H3K36me2	16-0319	50 µg
H3K36me3	16-0320	50 µg
H3K79me1	16-0367	50 µg
H3K79me2	16-0368	50 µg
H3K79me3	16-0369	50 µg
H4K20me1	16-0331	50 µg
H4K20me2	16-0332	50 µg
H4K20me3	16-0333	50 µg



dNucs

Other PTMs

H2AUb*	16-0020	50 µg
H2AK119ub <i>Coming Soon</i>	16-0363	50 µg
H2BK120ub	16-0370	50 µg
H3R2,8,17cit	16-0362	50 µg
H3S10ph	16-0345	50 µg

* Enzymatically-modified; contains ubiquitination at H2AK13/15 and H2AK119.



dNucs

Histone Arginine Methylation

H2AR3me1	16-0359	50 µg
H2AR3me2a	16-0360	50 µg
H2AR3me2s	16-0361	50 µg
H3R2me1	16-0340	50 µg
H3R2me2a	16-0341	50 µg
H3R2me2s	16-0355	50 µg
H4R3me1	16-0356	50 µg
H4R3me2a	16-0357	50 µg
H4R3me2s	16-0358	50 µg



dNucs

Histone Acylation

H3K4ac	16-0342	50 µg
H3K9ac	16-0314	50 µg
H3K9bu	16-0371	50 µg
H3K9cr <i>Coming Soon</i>	16-0351	50 µg
H3K14ac	16-0343	50 µg
H3K18ac	16-0372	50 µg
H3K18bu	16-0373	50 µg
H3K18cr	16-0337	50 µg
H3K23ac	16-0364	50 µg
H3K27ac	16-0365	50 µg
H4K5ac	16-0352	50 µg
H4K8ac	16-0353	50 µg
H4K12ac	16-0312	50 µg
H4K16ac <i>Coming Soon</i>	16-0354	50 µg
H3K4,9,14,18ac	16-0336	50 µg
H3K4me3, K9,14,18ac	16-0335	50 µg
H4K5,8,12,16ac	16-0313	50 µg

* Non-biotinylated available upon request

Other Recombinant Nucleosomes (Biotinylated)*



oncoNucs

AA Substitutions Implicated in Cancer

H3.3K4M	16-0349	50 µg
H3.3K9M	16-0350	50 µg
H3.3K27M	16-0323	50 µg
H3.3G34R <i>Coming Soon</i>	16-0346	50 µg
H3.3G34V	16-0347	50 µg
H3.3G34W	16-0348	50 µg
H3.3K36M	16-0344	50 µg



vNucs

Histone Variants

H2AX	16-0013	50 µg
H2AZ.1	16-0014	50 µg
H2AZ.2	16-0015	50 µg
H3.3	16-0011	50 µg
H3.3, non-biotinylated	16-0012	100 µg
H2AXS139phos	16-0366	50 µg



methyl DNA Nucs

Mononucleosomes, Hemi-methylated, Biotinylated	16-2003	50 µg
Mononucleosomes, Hemi-methylated	16-2103	50 µg
Mononucleosomes, 187x601 DNA, Biotinylated	16-2004	50 µg
Mononucleosomes, 187x601 DNA	16-2104	50 µg



rNucs

Human Recombinant, No PTMs

Mononucleosomes, biotinylated	16-0006	50 µg
Mononucleosomes, non-biotinylated	16-0009	100 µg

* Non-biotinylated available upon request

Custom nucleosome synthesis available.

Please email us at info@epicypher.com for additional information and pricing or complete a [Request for Quote](#) form on our website.

Functionalized Nucleosome Substrates for Drug Discovery and Chromatin Research

Recombinant Nucleosome Remodeling Substrates



EpiDyne® FRET Nucleosome Remodeling Assay Substrate **16-4201** 50 µg

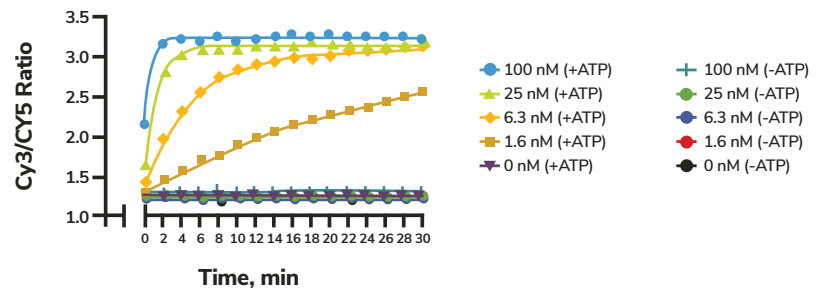
EpiDyne® Nucleosome Remodeling Assay Substrate ST601-GATC1 **16-4101** 50 µg

EpiDyne® Remodeling Assay Substrate DNA ST601-GATC0 **18-4100** 50 µg

EpiDyne® Remodeling Assay Substrate DNA ST601-GATC1 **18-4101** 50 µg

ATP-dependent nucleosome remodeling enzyme complex

ACF Remodeling Enzyme Complex **15-1013** 100 reactions

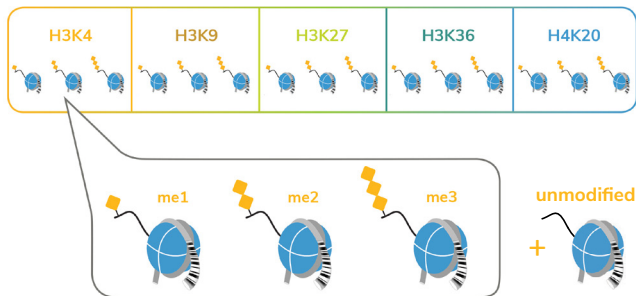


ACF Chromatin Remodeling Assay:

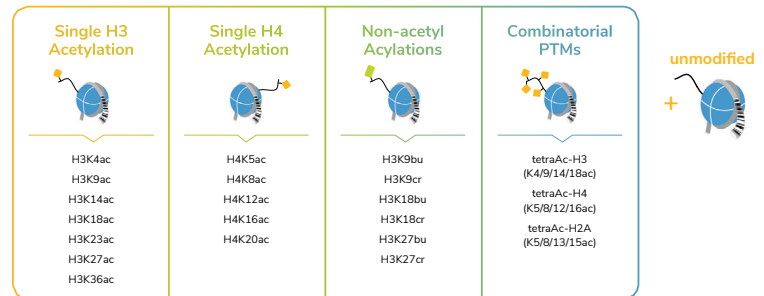
Assays were performed using EpiDyne®-FRET nucleosome remodeling substrate. EpiDyne®-FRET was incubated with increasing amounts of ACF Remodeling Enzyme Complex (as indicated), with or without ATP. ACF remodeling activity was measured by changes in FRET (Cy3/Cy5 ratio).

SNAP-CHIP® spike-in panels

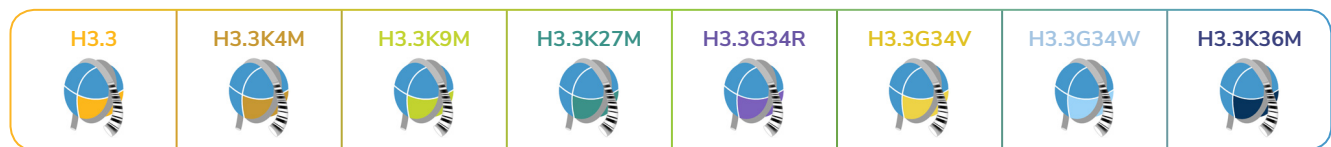
K-MetStat Panel (Cat. No. 19-1001)



K-AcylStat Panel (Cat. No. 19-3001)



Oncostat Panel (Cat. No. 19-2001)



ORDERING INFO

Let's discuss your project
sales@epicypher.com

For more information
epicypher.com/nucleosomes

To place an order
sales@epicypher.com

Coming Soon
R-MetStat panel

EpiCypher.com
855.374.2461
info@epicypher.com

