

dCypher® R-MetStat Nucleosome Panel

Catalog No. 16-9004
Lot No. 19070001
Pack Size 96 well plate



EpiCypher®

Product Description:

A 96-well plate containing 15 unique mononucleosomes assembled from recombinant human histones expressed in *E. coli* (two each of histones H2A, H2B, H3 and H4; accession numbers: H2A-P04908, H2B-O60814, H3.1-P68431/H3.2-Q71D13, H4-P62805). Single histone post-translational modifications (PTMs) are created using a proprietary synthetic method. The nucleosomes are wrapped by 147 base pairs DNA containing a central 601-positioning sequence (Lowary and Widom 1998 J Mol Biol. (1):19-42) with a 5' biotin-TEG group.

All nucleosomes in the panel are subjected to EpiCypher's rigorous quality control metrics, including: ESI-TOF mass spectrometry analysis of the modified proteins; SDS-PAGE to confirm octamer composition and purity; native PAGE to confirm nucleosome assembly and lack of free DNA; and Western blot analysis of the PTM. For the full list of nucleosomes in the panel, including individual catalog numbers of full-size (50 µg) products, see the associated excel sheet on the product page at www.epicypher.com.

	1	2	3	4	5	6	7	8	9	10	11	12
A	Unmodified	H3R8me2a			Unmodified	H3R8me2a			Unmodified	H3R8me2a		
B	H2AR3me1	H3R8me2s			H2AR3me1	H3R8me2s			H2AR3me1	H3R8me2s		
C	H2AR3me2a	H3R17me1			H2AR3me2a	H3R17me1			H2AR3me2a	H3R17me1		
D	H2AR3me2s	H3R17me2a			H2AR3me2s	H3R17me2a			H2AR3me2s	H3R17me2a		
E	H3R2me1	H4R3me1			H3R2me1	H4R3me1			H3R2me1	H4R3me1		
F	H3R2me2a	H4R3me2a			H3R2me2a	H4R3me2a			H3R2me2a	H4R3me2a		
G	H3R2me2s	H4R3me2s			H3R2me2s	H4R3me2s			H3R2me2s	H4R3me2s		
H	H3R8me1				H3R8me1				H3R8me1			
	Replicate #1		Blank		Replicate #2		Blank		Replicate #3		Blank	

dCypher® R-MetStat Nucleosome Panel plate layout. The R-MetStat panel includes an unmodified nucleosome control and 14 designer nucleosomes (dNucs: single PTMs, either monomethylation, dimethyl asymmetric, or dimethyl symmetric methylation at the indicated arginine residues).

Formulation:

Purified recombinant mononucleosomes individually stored in a 96 well plate (1.5 µg nucleosome at 1.5 µM per well) in 5 µL of nucleosome storage buffer (10 mM Tris-HCl pH 7.5, 1 mM EDTA, 25 mM NaCl, 2 mM DTT, & 20% glycerol). MW = ~200,000 Da

Storage and Stability:

Stable for six (6) months at -80°C from date of receipt. Avoid multiple freeze/thaws.

Application Notes:

The access to epigenetic diversity in a physiological nucleosome context enables broad end-user applications, including nucleosome binding studies (*e.g.* chromatin reader binding preferences), enzyme screening assays (*e.g.* identification of preferred substrates), and antibody specificity testing (*e.g.* for applications that require antibody specificity in a nucleosome, rather than histone peptide context). The biotin group on the DNA facilitates pull-down experiments. EpiCypher mononucleosomes do not contain free DNA which could alter assayed activities.

This product is for *in vitro* research use only and is not intended for use in humans or animals.

Frequently Asked Questions:

Q: *How many assay points of each nucleosome are supplied?*

A: This varies according to the sensitivity of the end application. For some context, the supplied format is sufficient to perform >35 AlphaScreen assays, >30 IP-qPCR experiments, and 3-5 immunoblotting experiments.

Q: *How do I store the plate? Can I dilute the nucleosomes?*

A: The nucleosomes are stable for 6 months at -80°C in the supplied format (1.5 µM). Freeze/thaw cycles should be limited (2-3 times). When diluted below 1.5 µM, a carrier protein (*e.g.* BSA) must be used to supplement the concentration, otherwise, stability is affected. For example, in our SNAP-ChIP product line, nucleosomes are diluted to 0.6 nM in 10 mM sodium cacodylate, pH 7.5, 100 mM NaCl, 1 mM EDTA, 50% glycerol (w/v), 1x Protease Inhibitor cocktail, 100 µg/mL BSA, 10 mM β-mercaptoethanol. In this formulation, the nucleosomes are stable at -20°C (where they do not freeze) for 6 months.

For applications that may not be compatible with the SNAP-ChIP buffer, EpiCypher routinely dilutes nucleosomes to 40 nM in assay buffer (50 mM Tris pH 7.5 + 0.01% Tween-20 + 0.01% BSA). Under these conditions the nucleosomes are stable for ~2 weeks at -20°C. *Note: extremely high amounts of salt (>650 mM) and ionic detergents (*e.g.* SDS) will disrupt nucleosome stability.

Q: *How can I use a few nucleosomes in the panel and avoid freeze/thaw of the rest?*

A: The nucleosomes are conveniently packaged in 8-well strip tubes and caps, so each column can be conveniently separated from the rest of the plate. For example, if the user desires to use only one of the three replicates in the plate, the tubes in columns 1 and 2 can be removed without thawing the remainder of the plate.

This product is for *in vitro* research use only and is not intended for use in humans or animals.