

## Product Information

<b>Description:</b>	Nucleosomes prepared from mammalian HeLa cells (depleted from Histone H1) as a substrate for histone methyltransferase assays.
<b>Background:</b>	Nucleosomes are portions of DNA that are wrapped around histone proteins in the form of a histone octamer. The histone octamer consists of two copies of H2A, H2B, H3 and H4. Nucleosomes correspond to subunits of chromatin and allow the DNA to be compacted enough to fit into the cell's nucleus. The presence of nucleosomes at specific points of the genome determines the ability of regulatory proteins to access DNA.
<b>Species</b>	Human
<b>Concentration:</b>	1.01 mg/ml
<b>Expression System:</b>	HeLa
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	20 mM Tris-HCl, pH 7.5, 1 mM EDTA, 3 mM DTT, 10 mM Na butyrate, and 20% glycerol
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Applications:</b>	Useful as substrate for histone methyltransferase assays.

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining

