

## Recombinant Histone H3 dimethyl Lys79

**Catalog No:** 31221

**Source:** *Xenopus laevis*

**Quantity:** 50 µg

**Expressed In:** *E. coli*

**Contents:** Lyophilized from 50 µl sterile water containing 5 mM Beta-mercaptoethanol. Supplied as lyophilized powder.

### Background:

**Histone H3** is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 146 base pairs of DNA wrapped around an octamer of core histone proteins (two each of H2A, H2B, H3 and H4). Histone H1 is a linker histone, present at the interface between the nucleosome core and DNA entry/exit points.

### Protein Details:

Recombinant *Xenopus laevis* Histone H3 dimethyl Lys79 produced in *E. coli* and purified using FPLC. The purified protein was subsequently alkylated to specifically **dimethylate lysine 79** and repurified prior to lyophilization. The degree of methylation is > 99% and confirmed by high-resolution ESI-TOF. Protein concentration was determined by Bradford assay. ≥ 98% purity by SDS-PAGE.

### Protein Applications:

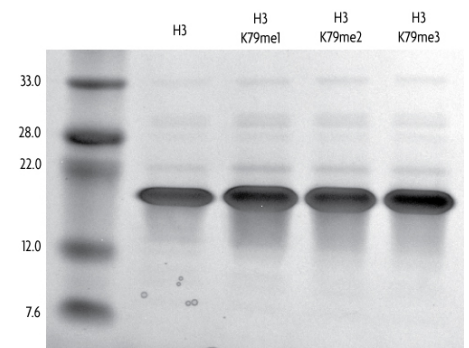
Each lot has been evaluated by SDS-PAGE and dot-blot analysis.

### Storage:

Lyophilized protein can be stored at -20°C. Once the protein is resuspended, we recommend making aliquots to avoid multiple freeze-thaw cycles and storing the protein at -80°C.

### Guarantee:

For research use only. Product is guaranteed stable for six months from date of receipt when stored properly.



SDS-PAGE analysis of 1.5 µg Recombinant Histone H3 (C110A) (lane 2), Recombinant Histone H3 monomethyl Lys79 (lane 3), Recombinant Histone H3 dimethyl Lys79 (lane 4), and Recombinant Histone H3 trimethyl Lys79 (lane 5). Molecular weight marker is in lane 1.