

## Histone H2AZ.1, Recombinant Human

<b>Catalog No</b>	15-0305	<b>Species</b>	Human
<b>Lot No</b>	15259001	<b>Source</b>	<i>E. coli</i>
<b>Pack Size</b>	100 µg	<b>Epitope Tag</b>	None
<b>Concentration</b>	N/A	<b>MW</b>	14 kDa

### DESCRIPTION

Recombinant human histone H2AZ.1 (H2A.Z; H2AFZ; P0C0S5) expressed in *E. coli* and purified by FPLC. Histone variants H2AZ.1 and H2AZ.2 (H2A.V; H2AF.V; Q71UI9) differ by three amino acids and are incorporated into a subset of nucleosomes in place of canonical histone H2A. H2AZ is implicated in diverse functions ranging from transcriptional regulation, chromosome transmission and DNA damage repair. H2AZ has a dedicated deposition machinery (the SWR-C ATPase) and is acetylated by the KAT5/TIP60 complex.

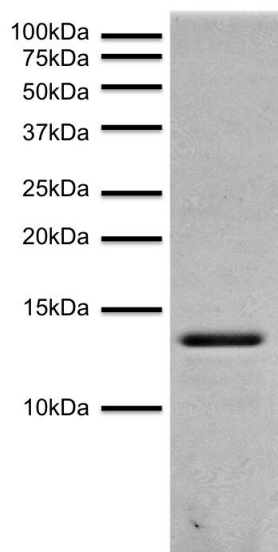
### TECHNICAL INFORMATION

<b>Storage</b>	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
<b>Formulation</b>	100 µg of lyophilized powder.

### APPLICATION NOTES

Recombinant histone H2AZ.1 is suitable for enzyme assays and nucleosome reconstitution. Reconstitute with distilled water or suitable buffer prior to usage.

### VALIDATION DATA



**FIGURE 1 Protein gel data.** Histone H2AZ.1 (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.