

Histone H2AZ.1, Recombinant Human

Catalog No. 15-0305
Lot No. 15259001
Pack Size 100 µg

Type Histone **Expressed In** *E. coli*
Mol. Wgt. 14 kDa **Epitope Tag** None

Product Description:

Recombinant human histone H2AZ.1 (H2A.Z; H2AFZ; POCOS5) 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.

Formulation:

100 µg of lyophilized powder.

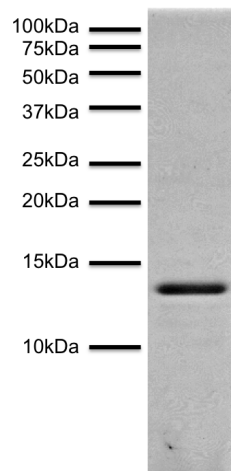
Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

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

References:



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.

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