

Histone H2AZ.2, Recombinant Human

Catalog No	15-0306	Species	Human
Lot No	16060002	Source	<i>E. coli</i>
Pack Size	100 µg	Epitope Tag	None
Concentration	N/A	MW	13.5 kDa

DESCRIPTION

Recombinant human histone H2AZ.2 (H2A.V; H2AF.V; Q71UI9) expressed in *E. coli* and purified by FPLC. Histone variants H2AZ.1 (H2A.Z; H2AFZ; P0C0S5) and H2AZ.2 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 KAT4/TIP60 complex.

TECHNICAL INFORMATION

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

APPLICATION NOTES

Recombinant histone H2AZ.2 is suitable for enzyme assays and nucleosome reconstitution. Reconstitute with distilled water prior to usage.

VALIDATION DATA

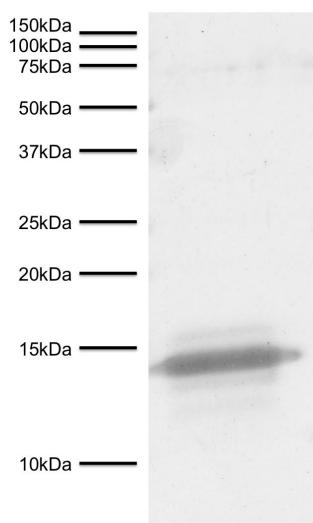


FIGURE 1 Protein gel data. Histone H2AZ.2 (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.