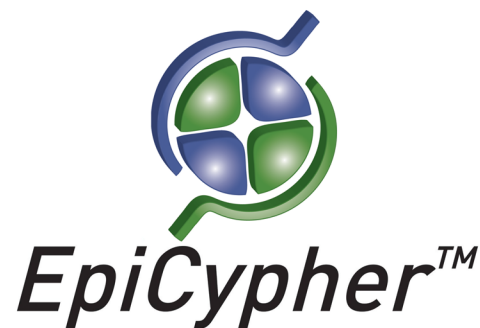


# Histone H2AZ.2, Recombinant Human

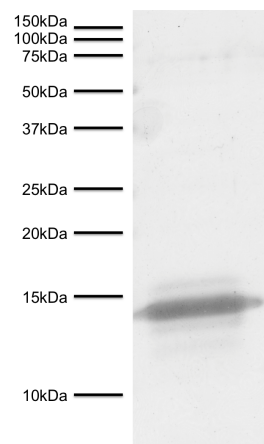
**Catalog No.** 15-0306  
**Lot No.** 16060002  
**Pack Size** 100 µg



**Type** Histone                      **Expressed In** *E. coli*  
**Mol. Wgt.** 13.5 kDa              **Epitope Tag** None

## Product 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; POC0S5) 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.



**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.

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

## References:

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