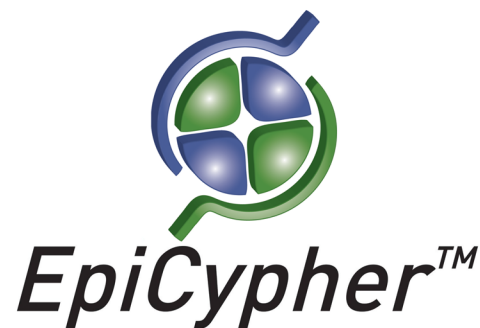


Histone H2AX, Recombinant Human

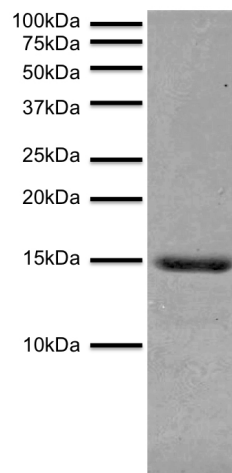
Catalog No. 15-0307
Lot No. 15259001
Pack Size 100 µg



Type Histone **Expressed In** *E. coli*
Mol. Wgt. 15 kDa **Epitope Tag** None

Product Description:

Recombinant human histone H2AX (H2AFX, accession P16104) expressed in *E. coli* and purified by FPLC. Histone H2AX is a histone variant that is incorporated into a subset of nucleosomes in place of canonical histone H2A. H2AX is phosphorylated on serine 139 by the ATM and ATR kinases in response to DNA double strand breakage. Phosphorylated H2AX serves as a platform to recruit a number of DNA damage repair proteins, including BRCA1, NBS1 and MDC1.



Protein Gel Data: Histone H2AX (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 H2AX is suitable for enzyme assays and nucleosome reconstitution. Reconstitute with distilled water or suitable buffer prior to usage.

References:

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