
H2AFX Protein

(Catalog No. E35001)

Description

H2AFX is one of several genes coding for histone H2A. H2AFX becomes phosphorylated on serine 139, then called gamma-H2AX, as a reaction on DNA Double-strand breaks (DSB). Human H2AFX partial ORF (AAH11694.1, 1 a.a. - 96 a.a.) recombinant protein with GST-tag at N-terminal.

Theoretical MW (kDa)

36.19

Preparation Method

In vitro wheat germ expression system

Size

50 µg

Storage Buffer

50 mM Tris-HCl, 10 mM reduced Glutathione, pH 8.0

Storage Instructions

Store at -80°C. Aliquot to avoid repeated freezing and thawing.

Ordering Information:

Products	Size	Cat. No.
H2AFX Protein	50 µg	E35001 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.