

Histone H3.X/Y antibody (mAb)

Catalog Nos: 61161, 61162

RRID: AB_2793533

Clone: 8H6-2111

Isotype: IgG2a

Application(s): ChIP, ICC, IF, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 17 kDa

Background: Histones H3.X and H3.Y are primate-specific variants of histone H3. H3.X and H3.Y are roughly 80% conserved compared to histone H3.3 The H3.Y protein is highly expressed in several primary cell lines, cancer cell lines and human brain. Expression of histone H3.Y is induced by stress. Exogenously expressed H3.X and H3.Y are efficiently incorporated into chromatin and exhibit dynamics similar to canonical histone H3. Knockdown of H3.Y results in slowed growth and altered expression of cell cycle related genes.

Immunogen: This H3.X/Y antibody was raised against a synthetic peptide corresponding to amino acids 9 - 20 of human H3.X.

Buffer: Purified rat IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.