

Histone H3K23pr antibody (pAb)

Catalog Nos: 61397, 61398

RRID: AB_2793619

Isotype: IgG

Application(s): DB, WB

Reactivity: Human, Wide Range Predicted

Volumes: 100 μ l, 10 μ l

Purification: None

Host: Rabbit

Molecular Weight: 17 kDa

Background: Propionylation of lysines is a recently identified post-translational modification and has been demonstrated to occur in proteins including histones and non-histone substrates, such as p53. **Propionylation of Lys23 of Histone H3** may be catalyzed by the histone acetyltransferase p300 and may be removed by histone deacetylase Sir2. H3K23 propionylation may play an important role in the epigenetic regulation of cell metabolism.

Immunogen: This antibody was raised against a synthetic peptide corresponding to propionyl-Lys23 human Histone H3.

Buffer: Rabbit serum containing 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.