
Histone H1.0 Monoclonal Antibody [RMC087A]

(Catalog # A68438)

Background

Histones H1 are necessary for the condensation of nucleosome chains into higher-order structures. The H1F0 histones are found in cells that are in terminal stages of differentiation or that have low rates of cell division.

Description

Histone H1.0 Monoclonal Antibody, clone RMC087A. Unconjugated. Raised in: Rabbit.

Concentration

[Inquire](#)

Formulation

Liquid. Rabbit IgG in phosphate buffered saline, pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Specificity

Human

Isotype

IgG

Uniprot ID

P07305

Purification

Affinity Purified

Immunogen

A synthesized peptide

Storage

Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Histone H1.0, Histone H1', Histone H1(0), Histone H1.0, N-terminally processed, H1F0, H1FV

Application

ELISA
