

# Histone H1

## Recombinant monoclonal antibody (HH1/1784R)

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. [UniProt]

This antibody is covered by our [Worry-Free Guarantee](#).

### Ordering Information

[Order Online »](#)

ENZ-ABS375-0100	100µg
-----------------	-------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## Product Details

Application	Flow Cytometry, IF, IHC (PS)
Clone	HH1/1784R
Formulation	Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
Host	Rabbit
Immunogen	Recombinant full length human Histone H1.
Isotype	IgGκ
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (0.5-1 µg/10 <sup>6</sup> cells in 0.1 mL)Immunofluorescence (0.5-1 µg/mL)Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 µg/mL for 30 min at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P07305
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>