

---

## Histone H4R3 Monomethyl (H4R3me1) Polyclonal Antibody

(Catalog #A-3717)

---

### Background

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. This structure consists of approximately 146 bp of DNA wrapped around a nucleosome, an octamer composed of pairs of each of the four core histones (H2A, H2B, H3, and H4). The chromatin fiber is further compacted through the interaction of a linker histone, H1, with the DNA between the nucleosomes to form higher order chromatin structures. This gene is intronless and encodes a member of the histone H4 family. Transcripts from this gene lack polyA tails; instead, they contain a palindromic termination element. This gene is found in a histone cluster on chromosome 1. This gene is one of four histone genes in the cluster that are duplicated; this record represents the centromeric copy.

### Description

Histone H4R3 Monomethyl (H4R3me1) Polyclonal Antibody, unconjugated. Raised in: Rabbit

### Formulation

PBS, pH 7.4, containing 0.02% sodium azide as preservative and 50% glycerol.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

P62805

### Purification

Affinity Purified

### Immunogen

A synthetic peptide

### Storage

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

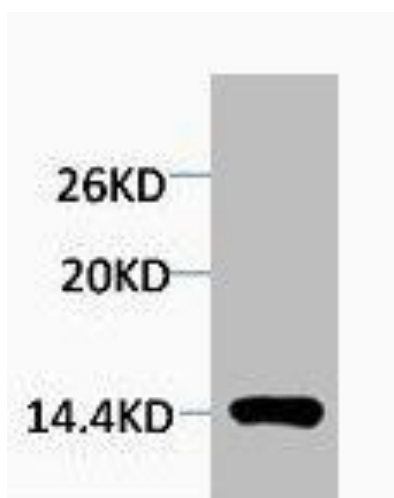
### Alternative Names

H4R3me1, H4R3m1, H4, H4/n, H4F2, H4FN, FO108, HIST2H4, H4 Arginine 3 me1

### Applications

WB; Recommended dilution: WB: 1:1000-2000

---



Western blot analysis of extracts from Hela cells, 1:2000