

Product Information

Description

Histone-H2A-(107-122)-Ac-OH is a peptide with the sequence Ac-Gly-Val-Leu-Pro-Asn-Ile-Gln-Ala-Val-Leu-Leu-Pro-Lys-Lys-Thr-Glu-OH, MW= 1762.1. Histone H2A is one of the five main histone proteins involved in the structure of chromatin in eukaryotic cells.

Target:

Cancer Biology

Pathway

Cancer Biology Pathway

Molecular Formula:

 $C_{81}H_{140}N_{20}O_{23}$

Supplied As:

Solid

Solubility:

≥176.2 mg/ml in DMSO; insoluble in EtOH; ≥23.3 mg/ml in H₂O. For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while. Long-term storage in solution is not recommended. Not stable in diluted solutions.

Stability:

At least 6 months at -20°C. Avoid freeze/thaw cycles.

Storage:

-20°C

MW:

1726.1 Da

Purity:

≥97%

Chemical Structure

