
SUV39H1 Protein

(Catalog No. E24037)

Description

Human SUV39H1 (GenBank Accession No. NM_003173), amino acids 82-412, with N-terminal GST tag.

Molecular Weight (kDa)

64 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCL, 3 mM DTT, 20 mM Glutathione and 10% Glycerol.

Source

Expressed in E.coli expression system

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity > 60%

Application Notes

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Ordering Information:

Products	Size	Cat. No.
SUV39H1 Protein	50 µg	E24037 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.