
SUV39H2 Protein

(Catalog No. E24038)

Description

Human SUV39H2 (GenBank Accession No. NM_024670), amino acids 26-350, with N-terminal GST tag

Molecular Weight (kDa)

63 kDa

Formulation

40mM Tris pH 8.0, 110mM NaCl, 2.2mM KCl, 16mM Glutathione 20% Glycerol, and 3mM DTT.

Source

Expressed in E.coli expression system

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

SDS-PAGE purified

Application Notes

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

Ordering Information:

| Products | Size | Cat. No. |
|-----------------|-------|-----------|
| SUV39H2 Protein | 50 µg | E24038 -1 |

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