
SET7 Protein

(Catalog No. E23000)

Description

Human SET7/9 recombinant histone methyltransferase with 366 amino acids.

Molecular Weight (kDa)

40.7KDa

Formulation

50 mM Tris-HCl buffer(pH7.5) 0.2M NaCl, 5mM DTT, 20%glycerol.

Source

Expressed in E.Coli expression system

Size

50 µg

Storage Temp & Conditions

Store at -20 C or -80 C for long term

Purification Notes

Purity: > 95%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
SET7 Protein	50 µg	E23000 -1

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