
SET8 Protein

(Catalog No. E24035)

Description

Human SET8 (GenBank Accession No. NM_020382), (amino acids 195-352) with N-terminal GST tag.

Molecular Weight (kDa)

44 kDa

Formulation

5 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 3 mM DTT, 18 mM glutathione and 10% glycerol.

Source

Expressed in an 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.
SET8 Protein	50 µg	E24035 -1

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