
G9a Protein

(Catalog No. E24032)

Description

Human G9a (euchromatic histone-lysine N-methyltransferase 2) (GenBank Accession No. NM_006709) (amino-acids 785-1210), with N-terminal GST tag.

Molecular Weight (kDa)

74 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM glutathione, 20% glycerol and 3mM DTT.

Source

Expressed in an E.Coli expression system.

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity>50%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
G9a Protein	50 µg	E24032 -1

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