
DOT1L Protein

(Catalog No. E24029)

Description

Human DOT1L (GenBank Accession No.NM_032482), amino-acids 2-416 with N-terminal GST tag, MW = 73 kDa, expressed in E.Coli expression system.

Molecular Weight (kDa)

73 kDa

Formulation

45 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 E.Coli expression system

Size

50 µg

Storage Temp & Conditions

>12 months at -80°C

Purification Notes

Purity >70%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
DOT1L Protein	50 µg	E24029 -1

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