
DNMT1 Protein

(Catalog No. E15000)

Description

Human DNMT1 (GenBank Accession No. NM_001130823), (a-a 2-1632), with N-terminal GST tag.

Molecular Weight (kDa)

211 kDa

Formulation

40mM Tris pH 8.0, 110mM NaCl, 2.2mM KCl, 0.04% Tween, 20% glycerol, and 3mM DTT.

Source

Expressed in Sf9 cells via a baculovirus expression system.

Size

10 µg

Storage Temp & Conditions

>6 months at -80°C. Avoid freeze/thaw cycles

Purification Notes

Glutathione Sepharose.

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
DNMT1 Protein	10 µg	E15000-1

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