
HDAC4 Protein

(Catalog No. E24007)

Description

Human HDAC4 (GenBank Accession No. NM_006037), amino acids 627- 1085 with N-terminal GST tag.

Molecular Weight (kDa)

75.2 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 18 mM glutathione, 10% glycerol.

Source

Expressed in baculovirus expression system.

Size

10 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >65%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
HDAC4 Protein	10 µg	E24007-1

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