
HDAC9 Protein

(Catalog No. E24012)

Description

Human HDAC9 (GenBank Accession No. NM_178423), amino acids 604- 1066 with C-terminal His tag

Molecular Weight (kDa)

50.7 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 250 mM imidazole 10% glycerol

Source

Expressed in baculovirus expression system

Size

10 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >80%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
HDAC9 Protein	10 µg	E24012-1

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