
HDAC1 Protein

(Catalog No. E24002)

Description

Human HDAC1 (GenBank Accession No. NM_004964), full length with C-terminal His-tag and C-terminal FLAG-tag.

Molecular Weight (kDa)

56 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol.

Source

Expressed in baculovirus expression system

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >60%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
HDAC1 Protein	50 µg	E24002-1

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