
HDAC2 Protein

(Catalog No. E24005)

Description

Human HDAC2 (GenBank Accession No. NM_001527), full length with C- terminal His-tag.

Molecular Weight (kDa)

56 kDa

Formulation

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

Source

Expressed in baculovirus expression system

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >94%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
HDAC2 Protein	50 µg	E24005-1

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