
HDAC5 Protein

(Catalog No. E24008)

Description

Human HDAC5 catalytic domain, variant 1 (GenBank Accession No.NM_005474), amino acid 657-1123 with C-terminal His-tag.

Molecular Weight (kDa)

51 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 225 imidazole, and 10% glycerol.

Source

Expressed in baculovirus expression system

Size

10 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >90%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
HDAC5 Protein	10 µg	E24008-1

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