
HDAC6 Protein

(Catalog No. E24009)

Description

Human HDAC6 (GenBank Accession No. BC069243), full length with N-terminal GST tag.

Molecular Weight (kDa)

159 kDa

Formulation

40 mM Tris pH 8.0, 346 mM NaCl, 20% glycerol, 0.8 mM DTT.

Source

Expressed in baculovirus expression system

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Affinity chromatography and gel filtration

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
HDAC6 Protein	50 µg	E24009-1

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