
HDAC10 Protein

(Catalog No. E24003)

Description

Human HDAC10 (a.a. 1-481) (GenBank Accession No. NM_032019) with N-terminal GST tag, MW= 79 kDa, expressed in baculovirus expression system.

Molecular Weight (kDa)

79 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 20 mM glutathione, and 10% glycerol.

Source

Expressed in baculovirus expression system

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Glutathione Sepharose

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
HDAC10 Protein	50 µg	E24003-1

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