
HDAC8 Protein

(Catalog No. E24011)

Description

Human HDAC8 (GenBank Accession No. NM_018486), full length with C-terminal His tag.

Molecular Weight (kDa)

46.4 kDa

Formulation

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

Source

Expressed in a baculovirus expression system

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >92%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
HDAC8 Protein	50 µg	E24011-1

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