
HDAC11 Protein

(Catalog No. E24004)

Description

Human HDAC11 (full length) (GenBank Accession No. NM_024827) with N-terminal GST tag.

Molecular Weight (kDa)

66 kDa

Formulation

40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM glutathione, and 20% 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.
HDAC11 Protein	50 µg	E24004-1

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