
HDAC3 Protein

(Catalog No. E24006)

Description

Complex of human HDAC3 (GenBank Accession No. NM_003883), full length with C-terminal His tag, and human NCOR2 (amino acid 395-489) (GenBank Accession No. NM_006312), N-terminal GST tag.

Molecular Weight (kDa)

HDAC3: 49.7 kDa

NCOR2: 37.6 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole and 20% glycerol

Source

Co-expressed in baculovirus expression system.

Size

50 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity: >91%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
HDAC3 Protein	50 µg	E24006-1

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