
HDAC10 Protein, FLAG-Tag Recombinant

(Catalog # E24040)

Description

Human HDAC10, also known as Histone Deacetylase 10, Polyamine Deacetylase HDAC10, and HD10, GenBank Accession No. NM_032019, a.a. 2-631 with N-terminal FLAG tag, expressed in a Baculovirus-infected Sf9 cell expression system.

Molecular Weight (kDa)

67 kDa

Formulation

40 mM Tris-HCl, pH 8.0. 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% Glycerol and 90 µg/ml FLAG peptide.

Source

Human species expressed in an Sf9 host species/expression system

Size

50 µg

Tag(s)

N-terminal FLAG-tag

Storage Temp & Conditions

Store at -80°C for up to 6 months. Avoid freeze/thaw cycles.

Purity

≥62%

Application Notes

Useful for investigating histone deacetylation, gene expression regulation, chromatin structure, and related biological processes.
