
UTX Protein

(Catalog No. E24028)

Description

Human UTX, also known as KDM6A (GenBank Accession No. NM_021140) (amino-acids 1-1401) with C-terminal FLAG-tag, MW = 155 kDa, expressed in Sf9 cells via a baculovirus expression system.

Molecular Weight (kDa)

155 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 90µg/ml, 10% glycerol, and 3 mM DTT.

Source

Expressed in Sf9 cells via a baculovirus expression system

Size

20 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity>75%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
UTX Protein	20 µg	E24028 -1

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