

Data Sheet

UTX (KDM6A), FLAG-tag

Human, Recombinant, N-terminal FLAG-tag

Catalog #: 50119

Lot #: 150806-3 **Conc.:** 0.23 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 80 µg/ml flag peptide and 0.2mM TCEP

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

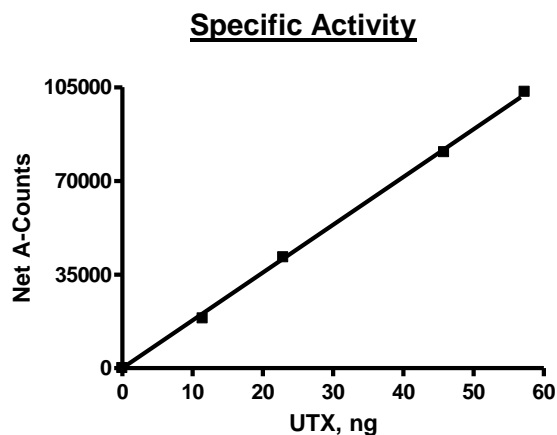
- Greenfield A, et al. (1998). *Hum. Mol. Genet.* **7** (4): 737–42.
- Agger K, et al. (2007). *Nature* **449** (7163): 731–4.

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

Specific Activity: ≥ 0.096 pmol/min/µg
Assay Conditions: 10 µl reaction mix containing assay buffer with 50 mM HEPES (pH 7.4), 1mM TCEP, 100 µM alpha-ketoglutarate, 25 µM iron, 0.1mM ascorbic acid, 0.02% Tween-20, 0.5 µM biotinylated peptide substrate, and UTX (10 – 50ng) added to the wells. Add antibody against dimethylated K27 peptide. Incubate for 30 min, then Streptavidin-conjugated secondary antibody followed by Alpha Screening detection.

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

Quality Assurance

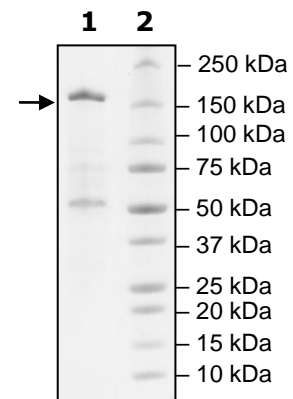


4-20% SDS-PAGE
Coomassie staining

Lane 1:
 UTX
 (1.2ug loaded)

Lane 2:
 Protein Marker

MW: 155 kDa
Purity: $\geq 58\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com