

KDM4A FL (StreptII) (Lysine-Specific Demethylase 4A, Full-length)

CATALOG NO.: PDM-21-437

LOT NO.:

DESCRIPTION: Full-length human recombinant KDM4A (residues 2-1064; Genbank Accession # NM_014663.2; MW = 122.9 kDa) expressed with an N-terminal StreptII tag in *sf9* insect cells.

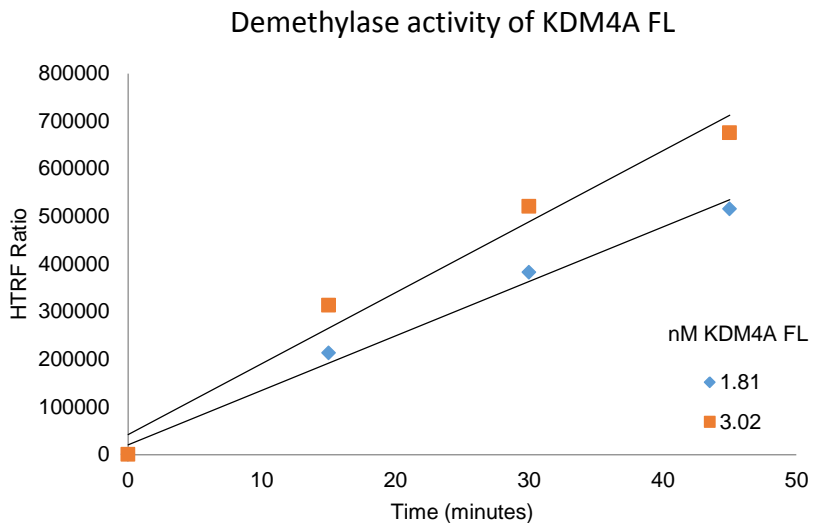
PURITY: >90% by SDS-PAGE

SUPPLIED AS: ___ µg/µL in 50 mM Tris HCl, pH 8, 300 mM NaCl, 5% glycerol, 0.5mM TCEP as determined by OD₂₈₀.

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 µg of RBC KDM4A FL (StreptII). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, 20, 15, & 10 kDa.



KDM4A FL Demethylase Activity Assay. B-H3(21-44)K36Me3 peptide demethylation was monitored using Cisbio HTRF detection (cat # 61KD2KAD). The 10 µL reaction contained 12.5 µM Fe(II), 20 µM α-KG, 100 µM ascorbate and 60 nM peptide and variable concentration of KDM4A FL. Reactions were quenched at specified time by the addition of detection reagent. Fluorescence emission (665 and 615 nm) was read using Envision reader (Perkin Elmer) following 3h incubation.

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.