

KDM5C (Strep) (Lysine-Specific Demethylase 5C)

CATALOG NO.: PDM-11-436

LOT NO.:

DESCRIPTION: Human recombinant KDM5C (residues 2-1560; Genbank Accession # NM_004187; MW = 178.6kDa) expressed with an N-terminal StrepII tag in *Sf9* insect cells.

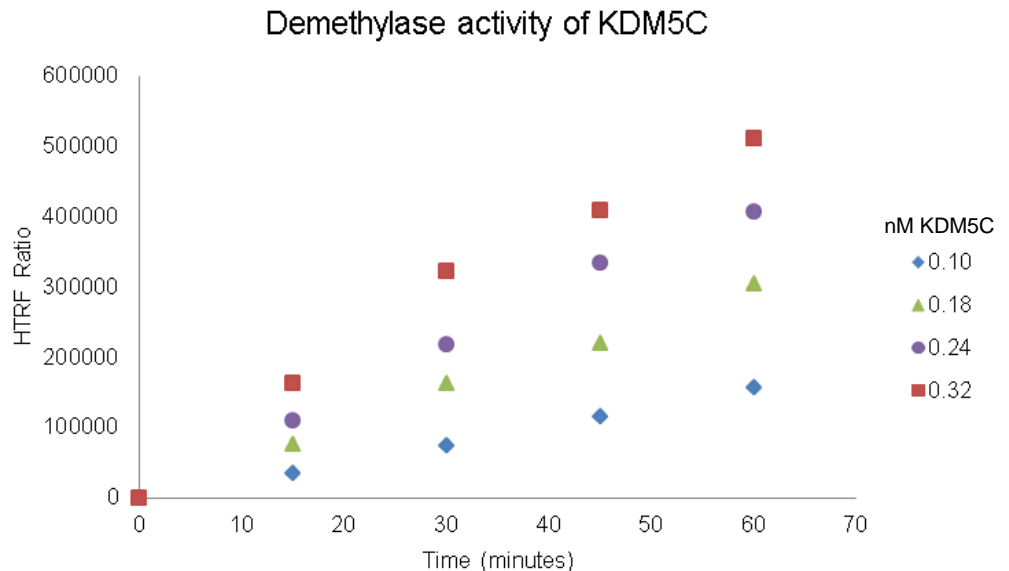
PURITY: >90% by SDS-PAGE

SUPPLIED AS: ___ $\mu\text{g}/\mu\text{L}$ in 50 mM Tris HCl, pH 8.0, 300 mM NaCl, 5% glycerol, 0.5 mM TCEP as determined by OD_{280} .

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 KDM5C (StrepII). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, & 10 kDa.



KDM5C Demethylase Activity Assay. B-H3(1-21)K4Me3 peptide demethylation was monitored using Cisbio HTRF detection (cat # 61KA2KAE). The 10 μL reaction contained 10 μM Fe(II), 15 μM α -KG, 100 μM ascorbate and 10 nM peptide and variable concentration of KDM5C. 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.