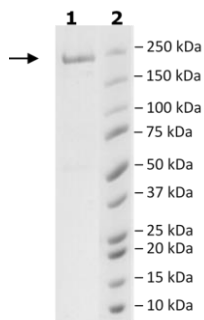


Product Information

Construct:	JMJD1B (KDM3B) (FLAG-Avi-2-1761(end))
Concentration:	0.72 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 80 µg/mL FLAG peptide
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_016604
MW:	195 kDa
Purity:	80%
Specific Activity:	0.02 pmol/min/µg
Assay Conditions:	Assay was performed according to the JMJD1B Homogeneous Assay Kit (BPS Bioscience #79839) with JMJD1B enzyme titrated from 400 ng/rxn – 0 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

