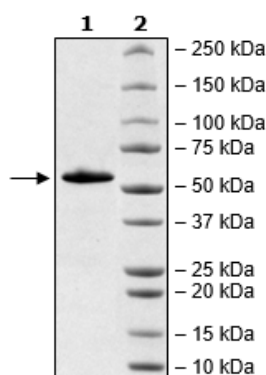


Product Information

Construct:	FTO (His-32-505(end))
Concentration:	0.56 mg/ml
Species:	Human
Formulated In:	40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 3 mM DTT
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001080432
MW:	56 kDa
Purity:	≥90%
Specific Activity:	0.003 pmole/min/μg
Assay Conditions:	Assay done using ELISA technology. 50 μl reaction mix containing HEPES-based buffer and 0-2000 ng/rxn FTO is added to the wells coated with the m6A substrate. Incubate overnight at room temperature. Add antibody against N6 methyladenosine, incubate 1 hr. Then, add secondary HRP-labeled antibody and incubate 30 min. Finally, add ELISA ECL substrates and read luminescence.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



FTO Activity

