

Data Sheet

FTO, His-Tag

Human, Recombinant, N-Terminal His-Tag

Catalog #: 79480

Lot #: 180412-1

Conc.: 0.76 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 20% glycerol,
 180 mM Imidazole

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and remain stable for 6 months at -80°C .

References:

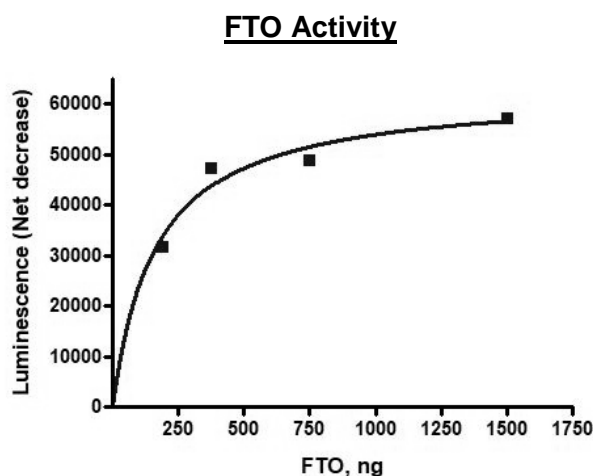
1. Frayling, T., *et al.*, *Science*. 2007; **316(5826)**: 889-894.
2. Dina, C., *et al.*, *Nature Genetics*. 2007; **39(6)**: 724-726.

Description: Human FTO, also known as Fat mass and obesity-associated protein, ALKBH9; BMIQ14; GDFD; KIAA1752, and alpha-ketoglutarate-dependent dioxygenase, GenBank Accession No. NM_001080432, a.a 32-505 (end), with N-terminal His-Tag and expressed in a Baculovirus infected Sf9 cell expression system. MW = 56 kDa.

Specific Activity: 0.0045 pmol/min/ μg
 Assay Conditions: Assay done using ELISA technology. 50 μl reaction mix containing HEPES-based buffer and 0-1500 ng FTO is added to the wells coated with the substrate. Incubate overnight at RT. Add antibody against N6-methyladenosine, incubate 1 hr. Then, add secondary HRP-labeled antibody and incubate 30 min. Finally, add HRP chemiluminescent substrates and read luminescence.

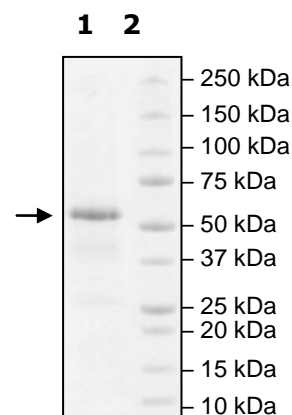
Applications: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 FTO
Lane 2:
 Protein Marker
MW: 56 kDa
Purity : $\geq 80\%$



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