

Data Sheet

FBXL10 (KDM2B, JHDM1B)

Human recombinant, with C-terminal FLAG-tag

Catalog #: 50120

Lot #: 140226

Conc: 0.21 mg/ml

Formulated in: 40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 ng/μL FLAG peptide, and 20% glycerol.

Stability: >6 months at -80°C

References:

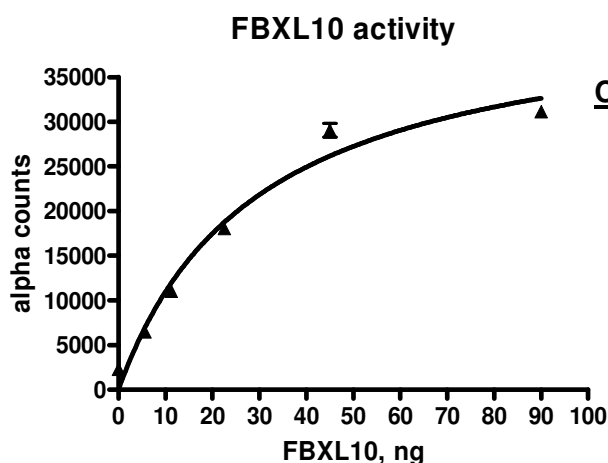
1. Tanaka, Y., *et al.* (2010). *EMBO J.* **29(9)**:1510-22.
2. Blackledge, N.P., *et al.* (2010). *Mol. Cell.* **38(2)**:179-90.
3. Frescas, D., *et al.* (2008). *Cell Cycle* **7(22)**:3539-47.

Description: Human F-Box Protein 10 (FBXL10, also known as KDM2B, JHDM1B, CXXC2, and Histone-H3 K36 demethylase 1B), GenBank Accession No. NM_032590, a.a. 1-650, with C-terminal FLAG-tag, MW = 75 kDa, expressed in Sf9 cells via a baculovirus expression system.

Assay Conditions: Enzyme was incubated with 0.4 μM substrate in HEPES-containing JMJ assay buffer SA and SA- (without KG) in a 10-μl reaction for 180 min at 30°C. 5 μl of Anti-Rabbit AlphaLISA® acceptor beads (1:250) and 5 μl of rabbit polyclonal antibody (1:200) were added and incubation continued for 30 min. Streptavidin Donor beads (1:125) were added (10 μl) and mixture was further incubated for 15 min at RT before A-counts were measured.

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

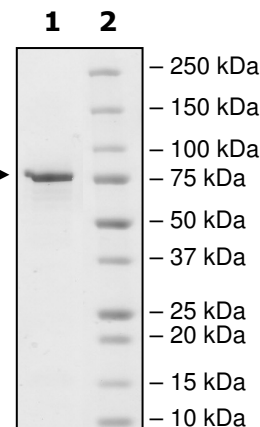
Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg FBXL10
Lane 2:
Protein Marker

MW: 75 kDa
Purity: ≥90%



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