

Data Sheet

ALKBH5, FLAG-Tag

Human, Recombinant, N-terminal FLAG-Tag

Catalog #: 100057

Lot #: 180723 **Conc.:** 0.56 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3mM DTT, 100ug/ml flag peptide

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Zheng, G. *et al.*, *Molecular Cell*. Jan; **46(1)**: 18-29 (2013).
2. Zhang, C. *et al.*, *PNAS*. **113(14)**: 47-56 (2016).

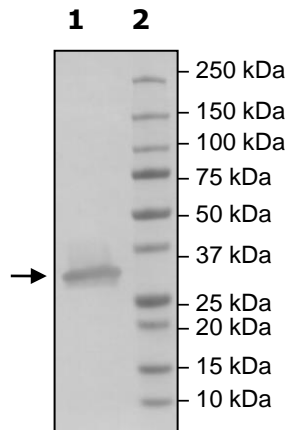
Description: Human ALKBH5, also known as AlkB Homolog 5, and Alkylated DNA repair protein alkB homolog 5, a.a. 66-292, with N-terminal FLAG-tag, GenBank Accession No. NM_017758, MW = 27 kDa, expressed in an Sf9 expression system.

Application: Useful as a positive control for Western blot studies.

Quality Assurance

**Western Blot
w/ FLAG-anti body**

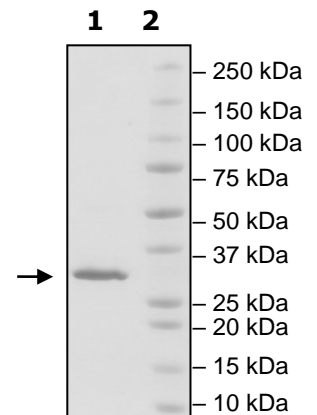
Lane 1:
ALKBH5
Lane 2:
Protein Marker



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

Lane 1:
ALKBH5
Lane 2:
Protein Marker

Purity: ≥90%
MW:
ALKBH5: 27 kDa



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