

Data Sheet

METTL3, FLAG-Tag

Human, Recombinant, N-terminal FLAG-tag

Catalog #: 100055

Lot #: 180626 **Conc.:** 0.077 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. For optimal
 storage, aliquot target into smaller
 quantities after centrifugation and store at
 recommended temperature.

References:

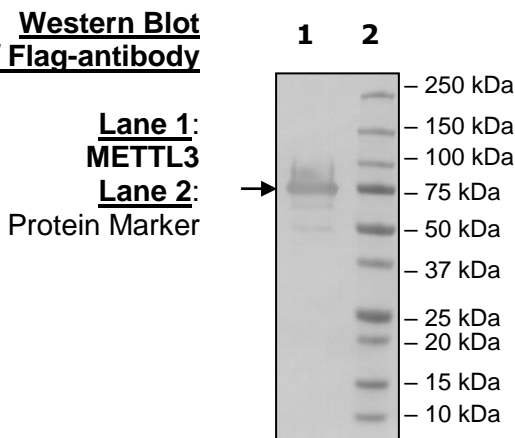
1. Liu J. *et al. Nat. Chem. Biol.* 2014. **10**: 93–95
2. Wang X. *et al. Nature.* 2016. **534**:575-578

Description: Human Methyltransferase-like 3 (METTL3), GenBank Accession No. NM_019852, 2-580 (end) with N-terminal FLAG-tag, expressed by baculovirus in Sf9 insect cells. MW = 66 kDa.

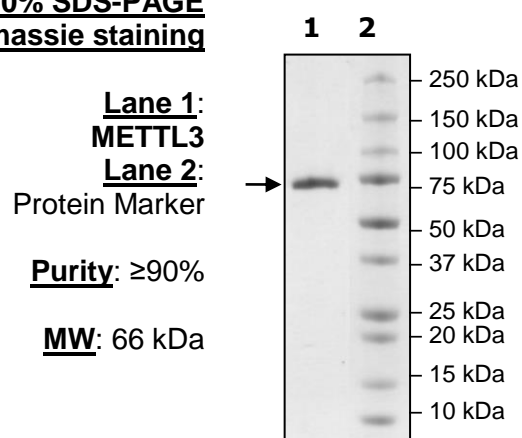
Application: Useful as a positive control for Western blot studies.

Quality Assurance

Western Blot
w/ Flag-antibody



4-20% SDS-PAGE
Coomassie staining



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