

## Data Sheet

### Lin28B, FLAG-Avi-Tag, Biotin-Labeled

Human, Recombinant, N-terminal FLAG-Avi-tag,  
Biotin Labeled

**Catalog #:** 79073

**Lot#:** 170711

**Conc.:** 0.13 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 100µg/ml FLAG  
peptide, 3mM DTT and 20% glycerol.

**Description:** Human LIN28B also known as  
Lin-28 Homolog B, with N-terminal FLAG-  
Avi-Tag and biotin labeled, GenBank  
Accession No. NM\_001004317, a.a. 2-250  
(end), expressed in a HEK293 cell  
expression system. MW = 30 kDa.

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

#### References:

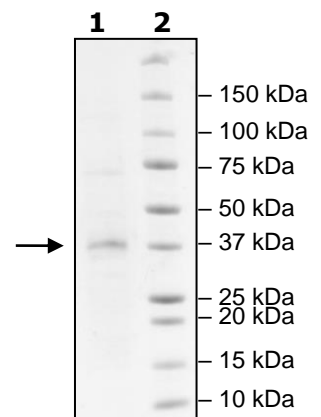
1. Piskounova, E. *et al.*, *Cell*. 2011; **5(23)**:  
1066-1079.
2. Tommiska, J. *et al.*, *BMC Research  
Notes*. 2011; **4(363)**.

---

### Quality Assurance

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

**Lane 1:**  
1µg Lin28B  
**Lane 2:**  
Protein Marker  
**MW:** 30 kDa  
**Purity:**  $\geq 50\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC USE.

To place your order, please contact us by Phone **1.858.829.3082**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)