

## Data Sheet

### Sestrin1

Human, recombinant, N-terminal FLAG-tag

Catalog #: 50284

Lot#: 100427

Conc.: 0.31 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 3 mM DTT, 90 µg/ml FLAG peptide and 10% glycerol.

**Stability:** >6 months at -80°C

#### References:

1. Lee, J.H., *et al.* (2010). *Science* **327**(5970): 1223-8.
2. Broderick, G., *et al.*, (2006). *Pharmacogenomics* 7(3): 407-419.
3. Budanov AV, Karin M. (2008). *Cell* **134** (3): 451-460.
4. Hay, N. (2008). *Cell Metab.* **8**(3):184-5.

#### Description:

Sestrin1, also known as p53-regulated protein PA26 (GenBank Accession No. NM\_014454), full length, a.a. 2-551(end), with N-terminal FLAG-tag, MW=65 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

#### Application:

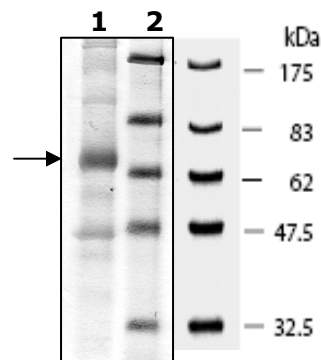
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance

#### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
4 µg Sestrin1  
**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 65 kDa  
**Purity:** >44%



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