

Data Sheet

Sestrin2

Human, recombinant, N-terminal FLAG-tag

Catalog #: 50286

Lot#: 100427

Conc.: 0.2 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:

Sestrin2, also known as hypoxia induced gene 95 (GenBank Accession No. NM_031459), full length, a.a. 2-480(end), with N-terminal FLAG-tag, MW=55 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

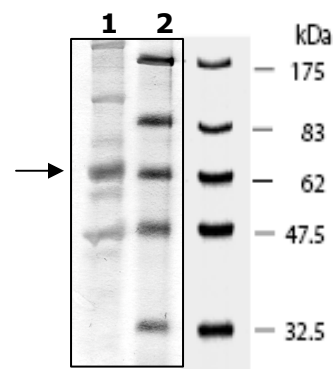
10% SDS-PAGE Coomassie staining

Lane 1:
3 µg Sestrin2

Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 55 kDa

Purity: >35%



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