

Data Sheet

SUZ12

Human, recombinant, His, Flag-tag

Catalog #: 50282

Lot#: 120912-C **Conc.:** 0.26 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.05% Tween-20, 20% glycerol, 1 mM DTT, and 100 µg/mL Flag peptide.

Stability: >6 months at -80 °C

References:

1. Cao, R, Zhang, Y. *Mol Cell*. **15(1)**: 57-67 (2004).
2. Brecqueville, M, *et al. Blood Cancer J*. **1(8)**:e33 (2011).
3. Li, H., *et al. Mol Cancer Res*. 2012 Sep 10. [Epub ahead of print]

Description:

Human SUZ12 (suppressor of zeste 12), Genbank Accession #NM_015355, (a.a. 2-739 (end)) with N-terminal His and Flag tags, MW = 85 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Application:

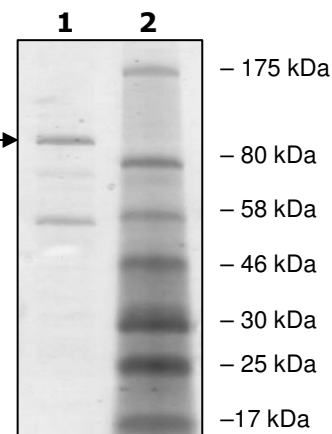
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% gradient SDS-PAGE Coomassie staining

Lane 1:
2 µg SUZ12
Lane 2:
Protein Marker

MW: 85 kDa
Purity: ≥65%



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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MHHHHHHDYKDDDDKAPQKHGGGGGGSGPSAGSGGGGFGGSAAVAAATASGG
KSGGGSCGGGGSYSASSSSSAAAAAGAAVLPVKKPKMEHVQADHELFLQAFEKPT
QIYRFLRTRNLIAPIFLHRTLTYMSHRNSRTNIKRKTFKVDDMLSKEKMKGEQESHSL
SAHLQLTFTGFFHKNDKPSNSENEQNSVTLEVLLVKVCHKKRKDVSCPIRQVPTGK
KQVPLNPDLNQTKPGNFPSLAVSSNEFEPNSHMKVSYLLFRVTRPGRREFNGMIN
GETNENIDVNEELPARRKRNREDGEKTFVAQMTVFDKNRRLQLLDGEYEVAMQEME
ECPISKKRATWETILDGKRLPPFETFSQGPTLQFTLRWTGETNDKSTAPIAKPLATRN
SESLHQENKPGSVKPTQTIAVKESLTTDLQTRKEKDPNENRQKLRIFYQFLYNNNTR
QQTEARDDLHCPWCTLNCRKLYSLLKHLKLCHSRFIFNYVYHPKGARIDVSINECYDG
SYAGNPQDIHRQPGFAFSRNGPVK RTPITHILVCRPKRTKASMSEFLESEGEVEQQR
TYSSGHNRLYFHSDTCLPLRPQEMEVDSEDEKDPEWLREKTITQIEEFSDVNEGEKEV
MKLWNLHVMKHGFIADNQMNHACMLFVENYGQKIIKKNLCRNFMHLVSMHDFNLISI
MSIDKAVTKLREMQQKLEKGESASPANEEITEEQNGTANGFSEINSKEKALETDSVSG
VSKQSKKQKL

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