

## Data Sheet

### ANGPT2

Human, recombinant, N-terminal FLAG-tag  
Catalog #: 91003  
Lot #: 130107                      Conc.: 0.06 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 100 ng/μl Flag peptide, and 20% glycerol.

**Stability:** >6 months at -80 °C

**References:**

1. Fiedler, U., *et al.*, *J. Biol Chem.* 2003 Jan 17;**278(3)**:1721-17.
2. Yuan, H.T., *et al.*, *Mol Cell Biol.* 2009 Apr;**29(8)**:2011-22.

**Description:**

Human angiopoietin 2 (ANGPT2), also known as ANG-2, GenBank Accession No. NM\_001147, a.a. 258-496 (end), with N-terminal FLAG-tag and mouse Cartilage Oligomeric Matrix Protein (COMP), GenBank Accession No. NM\_016685, a.a. 28-72, MW=33.4 kDa, expressed in a HEK293 cell expression system.

**Application:**

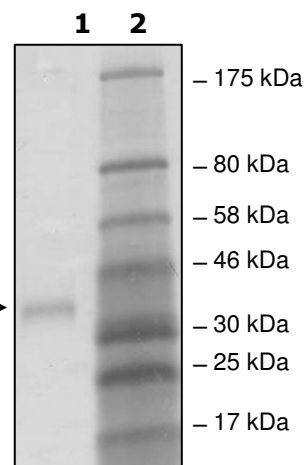
Useful for receptor binding studies, screening inhibitors, and selectivity profiling.

## Quality Assurance

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

**Lane 1:**  
1 μg ANG2  
**Lane 2:**  
Protein Marker

**MW:** 33.4 kDa  
**Purity:** ≥95%



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)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Sequence:**

DYKDDDDKDLAPQMLRELQETNAALQDVRELLRQQVKEITFLKNTVMECDACGLLTM  
MSTSNSAKDPTVAKEEQISFRDCAEVFKSGHTTNGIYTLTFPNSTEEIKAYCDMEAGG  
GGWTIIQRREDGSVDFQRTWKEYKVGFGNPSGEYWLGNFVSQLTNQQRYVLKIHKL  
DWEGNEAYSLYEHFYLSEELNYRIHLKGLTGTAGKISSISQPGNDFSTKGDNDKCI  
CKCSQMLTGGWWFDACGPSNLNGMYYPQRQNTNKFNGIKWYYWKSGYSLKATT  
MMIRPADF

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)