

Data Sheet

ANGPT1

Human, recombinant, N-terminal FLAG-tag
Catalog #: 91002
Lot #: 121113 Conc.: 0.23 mg/ml

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

Stability: >6 months at -80 °C

References:

1. Davis, S., *et al.*, *Cell*. 1996 Dec 27;**87**(7):1161-9.
2. Fiedler, U., *et al.*, *J. Biol Chem*. 2003 Jan 17;**278**(3):1721-17.

Description:

Human angiopoietin 1 (ANGPT1), also known as ANG-1, GenBank Accession No. NM_001146, a.a. 256-498 (end), with N-terminal FLAG-tag and mouse Cartilage Oligomeric Matrix Protein (COMP), GenBank Accession No. NM_016685, a.a. 28-72, MW=34 kDa, expressed in a HEK293 cell expression system.

Application:

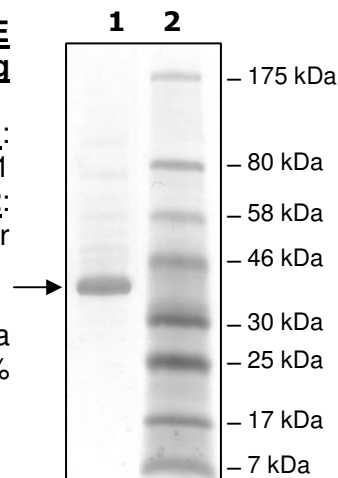
Useful for receptor binding studies, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE
Coomassie staining

Lane 1:
3 μg ANG1
Lane 2:
Protein Marker

MW: 34 kDa
Purity: ≥90%



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 W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

DYKDDDDKDLAPQMLRELQETNAALQDVRELLRQQVKEITFLKNTVMECDACGDTVHN
LVNLCTKEGVLLKGGKREEEKPFRCADVQAGFNKSGIYTIYINNMPKPKVFCNMD
VNGGGWTVIQHREDGSLDFQRGWKEYKMGFGNPSGEYWLGNEFIFAITSQRQYMLR
IELMDWEGNRAYSQYDRFHIGNEKQNYRLYLKGHTGTAGKQSSLILHGADFSTKDAD
NDNCMCKCALMLTGGWWFDACGPSNLNGMFYTAGQNHGKLNGIKWHYFKGPSYSL
RSTTMMIRPLDF

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