

# Angiopoietin-1 (human), (recombinant)

ANGPT1 is an angiogenic factor which intervenes in blood vessel maturation and takes part in endothelial development. ANGPT1 is a secreted ligand for Tie-2, a cell surface receptor tyrosine kinase expressed in endothelial and hemopoietic cells. The glycosylated ANGPT1 protein has a coiled-coil region in the amino terminus and a fibrinogen-like domain at the carboxy terminus.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-201-850-0020	20µg
------------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Store at +4°C if entire vial will be used within 2-4 weeks. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Ang-1, ANGPT1
<b>Appearance</b>	Colorless liquid.
<b>Concentration</b>	~1mg/ml
<b>Formulation</b>	Liquid. Sterile filtered, in 20mM TRIS-HCl buffer, pH 8.0, 1M urea containing 5% glycerol.
<b>MW</b>	~55.6kDa
<b>Purity</b>	≥85% (SDS-PAGE)
<b>Source</b>	Produced in <i>E. coli</i> . Human Ang-1 (angiopoietin-1) is a single, non-glycosylated polypeptide chain containing 480 amino acids (20-498)
<b>UniProt ID</b>	Q15389



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)