

COMP (rat):Angiopoietin-1 (human), (recombinant)

Soluble, stable and more potent Ang-1 variant.
Soluble, stable and more potent Ang-1 variant.

Citations: 12

[View Online »](#)

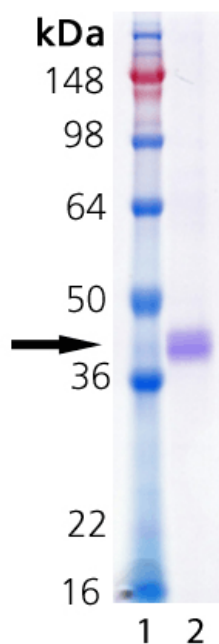
Ordering Information

[Order Online »](#)

ALX-201-314-C010	10µg
ALX-201-314-C050	50µg
ALX-201-314-C500	500µg

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis of ALX-201-314: Lane 1: MW Marker, Lane 2: 1µg of COMP (rat):Angiopoietin-1 (human), (recombinant) protein.

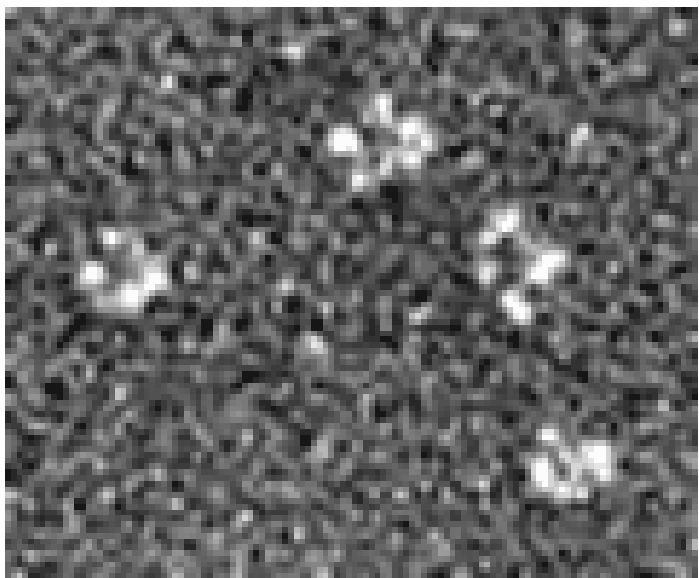


Figure 1: COMP (rat):Angiopoietin-1 (human) (rec.) (Prod. No. [ALX-201-314](#)) forms mainly pentameric structures. **Method:** Glycerol spraying/low angle rotary metal-shadowed specimens were imaged by transmission electron microscopy. COMP (rat):Angiopoietin-1 (human)



Structural scheme of ALX-201-314: Produced in CHO cells. The fibrinogen-like domain (aa 284-498) of human Ang-1 (angiopoietin-1) is fused at the N-terminus to the coiled-coil domain of rat COMP (cartilage oligomeric matrix protein) and to a FLAG®-tag.

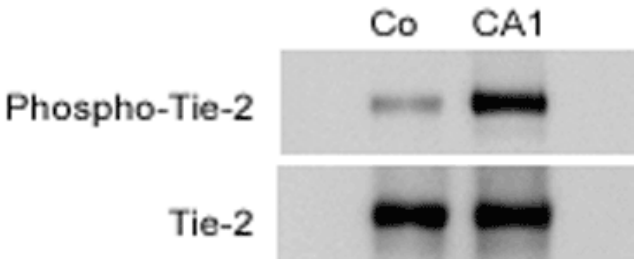


Figure 2: COMP (rat):Angiopoietin-1 (human) (rec.) (Prod. No. [ALX-201-314](#)) induces phosphorylation of Tie-2 in human umbilical vein endothelial cells. **Method:** Serum-starved human umbilical vein endothelial cells were treated with 200ng/ml of BSA (Co) or 200ng/ml of COMP:Angiopoietin-1 for 15 min. The phosphorylation of Tie-2 (phospho-Tie-2) was measured after immunoprecipitation. COMP (rat):Angiopoietin-1 (human)

Handling & Storage

Use/Stability	Working aliquots are stable for up to 3 months when stored at -80°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	COMP:Ang-1
Biological Activity	Induces phosphorylation of Tie-2 in primary cultured human and mouse vascular endothelial cells. More potent than native Ang-1.
Concentration	~0.2-0.4 mg/mL(Please refer to Certificate of Analysis for lot-specific concentration.)
Endotoxin Content	<0.1EU/μg protein (LAL test; BioWhittaker)
Formulation	Liquid. 0.2μm-filtered solution in PBS containing 0.01% CHAPS.
Purity	≥90% (SDS-PAGE)
Source	Produced in CHO cells. The fibrinogen-like domain (aa 284-498) of human Ang-1 (angiotensin-1) is fused at the N-terminus to the coiled-coil domain of rat COMP (cartilage oligomeric matrix protein) and to a FLAG®-tag.
Specificity	Binds to human Tie-2.

Technical Info / Product Notes FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID Q15389 (human Angiotensin-1), P35444 (rat COMP)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024