



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

Human Angiopoietin-2 (Ang2)

Catalog #: 91007
Lot #: 141103

Description: Recombinant human angiopoietin-2 (Ang2), also known as Angpt2, expressed in HEK293 cells. Human Ang2 is a 70 kDa secreted ligand that binds the receptor Tie2.

Source: Optimized DNA sequence encoding the extracellular domain of human Ang2, including a C-terminal His tag expressed in HEK293 cells.

Formulation: Recombinant human Ang2 is lyophilized from 0.2 μ m filtered 20 mM PBS solution, pH 7.2, 5% Trehalose.

Reconstitution: Briefly centrifuge the vial, followed by reconstitution in distilled water to a concentration of \geq 0.1 mg/ml. This solution can then be further diluted into other buffers containing carrier protein (0.1% BSA).

Storage: The lyophilized protein is stable for at least 1 year from date of receipt when stored at -20°C to -80°C. Upon reconstitution, store in working aliquots at 2 - 8°C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

Purity: >97%, as determined by SDS-PAGE and HPLC.

Endotoxin Level: Endotoxin level was found to be < 0.1 ng/ μ g (1EU/ μ g), using the LAL gel clot method.

Biologic Activity: The activity was tested by the ability of immobilized recombinant Human Tie2 to bind biotinylated recombinant human Ang2 in ELISA.

References:

1. *Circ. Res.* 2013; **113**: 1331-1344
2. *Science.* 2014; **343**: 1639-1646.

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