

Data Sheet

Rabbit Angiopoietin 2, His-tag

Rabbit, recombinant, C-terminal His-tag

Catalog #: 91010

Lot #: 150204-D1 **Conc.:** 0.54 mg/ml

Formulated in: 8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

Stability: At least 6 months at -80°C . *Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

Endotoxin Level <1 EU/ μg

References:

1. Mazziere, R., *et al.*, *Canc. Cell.* 2011; **19(4)**: 512-526.
2. Kim, I., *et al.*, *Oncogene.* 2000; **19(39)**: 4549-4552.

Description: Rabbit Angiopoietin 2 (Ang2), also known as ANGPT2, GenBank Accession No. XM_002720896.2, a.a. 19-496 with C-terminal His tag, MW=57.7 kDa, expressed in a HEK293 cell expression system. *Note:* this protein is expressed with a signal sequence (a.a. 1-18) that is removed after the protein is secreted. This protein runs at an apparent higher MW on SDS PAGE due to glycosylation.

Application: Useful for studies of protein function and protein-protein interaction.

Quality Assurance

4-20% SDS-PAGE
Coomassie staining

Lane 1:

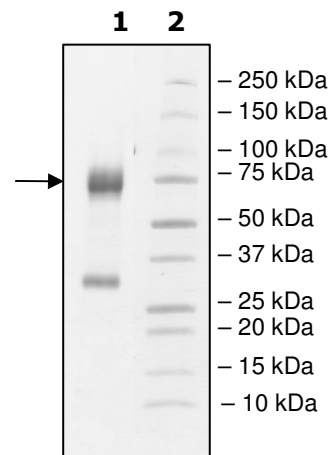
3 μg Ang2

Lane 2:

Protein Marker

MW: 57.7 kDa + glycans

Purity: $\geq 70\%$



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