

Data Sheet

Angiopoietin-like 3, FLAG-tag

Human recombinant with C-terminal FLAG-tag

Catalog #: 91009-2

Lot #: 141112

Amt: 50 µg

Formulated in: Lyophilized powder from 8 mM phosphate, pH 7.4, 100 mM NaCl, and 2.2 mM KCl

Stability: At least 12 months at -80°C as supplied.

Purity: ≥90% as determined by SDS-PAGE

Endotoxin Level: < 1 EU/µg

References:

1. Ando, Y., *et al.*, *J. Lipid Res.* **44**: 1216-1223.
2. Lee, E-Chiang, *et al.*, *J. Biol. Chem.* **284**: 13735-13745.

Description: Human angiopoietin-like 3 (ANGPTL3) also known as angiopoietin-related protein 3, or angiopoietin 5 (Ang5), GenBank Accession No. NM_014495.3, a.a. 1-460 with a C-terminal FLAG tag, expressed in a HEK293 cell expression system.

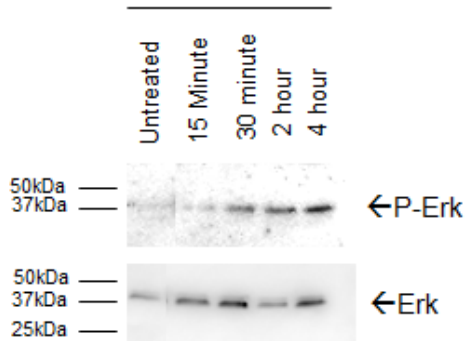
Biological Activity: Angiopoietin-like 3 activity was confirmed by treating THP-1 cells with 500 ng/ml of angiopoietin-like 3 for 15 min, 30 min, 2 hr, or 4 hr. Phosphorylated-Erk was measured by Western blot and compared to total Erk levels.

Application: Useful for *in vitro* cell culture assays or *in vivo* applications.

Quality Assurance

Biological Activity

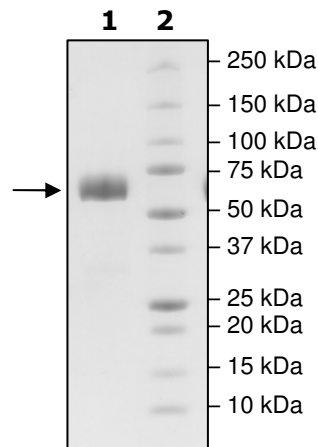
ANGPTL3 500 ng/ml



4-20% SDS-PAGE Coomassie staining

Lane 1:
3 µg ANGPTL3
Lane 2:
Protein Marker

MW: 52.8kDa
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