

## Data Sheet

### Angiopoietin-like 3, FLAG-tag

Human recombinant with C-terminal FLAG-tag

**Catalog #:** 91009-1

**Lot #:** 141112

**Amt:** 10 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.

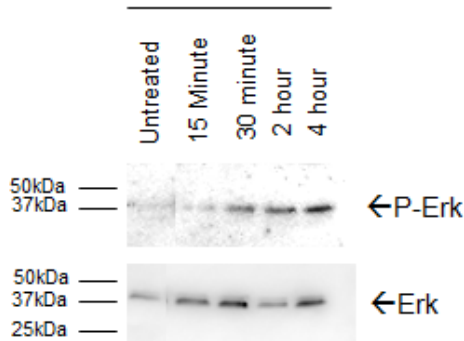
**Biological Activity:** Angiopoietin-like 3 activity was confirmed by treating THP-1 cells with 500 ng/ml of angiopoietin-like 3 for 15 min, 30 min, 2 hr, or 4 hr. Phosphorylated-Erk was measured by Western blot and compared to total Erk levels.

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

## Quality Assurance

### Biological Activity

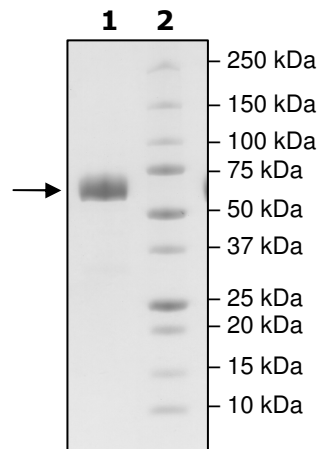
ANGPTL3 500 ng/ml



### 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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)