

## Data Sheet

### Nedd 4

Human, recombinant, N-terminal FLAG tag

**Catalog #:** 80404

Lot#: 120416-A      Conc.: 0.21 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 µg/ml Flag peptide, 20% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80°C

**References:**

1. Kotorashvili A. *et al.*, *J. Biol. Chem.* **284** (24), 16667-16678 (2009).
2. Plant P.J. *et al.*, *Biochem. J.* **419** (1), 57-63 (2009).

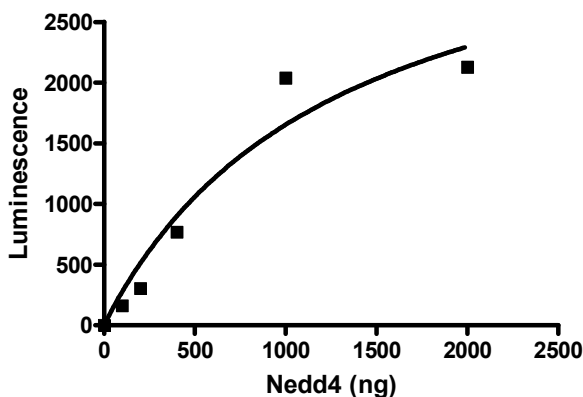
**Description:** Human Nedd 4 (GenBank Accession No. NM\_006154), amino-acids 150-end, with N-terminal FLAG tag, MW= 86 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Assay conditions:** 50 µl of reaction mix (4 mM ATP, 5 mM MgCl<sub>2</sub>, 1 mM DTT, E1 (0.5 µg), E2 (2.8 µg), Biotin-Ubiquitin (0.5 µg) and Nedd 4 (0-1.25 µg) in 50 mM Tris pH 7.4) is incubated for 30 min at 37°C. Transfer the reaction mixture to a streptavidin-coated plate and incubate for 30 min at 37°C. Add antibody against Nedd 4, incubate 1 hr. Then, add secondary HRP-labeled antibody and incubate 30 min. Finally, add HRP chemiluminescent substrates and read luminescence.

**Application:** Useful in conjunction with E1 and E2 for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance

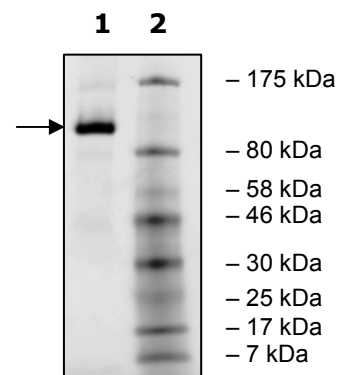
**Nedd 4 Specific Activity**



**10% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
 3.2 µg Nedd 4  
**Lane 2:**  
 Protein Marker  
 BioLabs (#P7708L)

**MW:** 86 kDa  
**Purity:** ≥92%



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)