

## Data Sheet

### SMURF 1, FLAG-Tag

Human, Recombinant, N-terminal FLAG tag  
**Catalog #:** 80402  
**Lot#:** 160314-G2      **Conc.:** 0.27 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Protein may be diluted to  $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored at  $-80^{\circ}\text{C}$ .

**References:**

1. Narimatsu M. *et al.*, *Cell* **137** (2), 295-307 (2009)
2. Fukunaga E. *et al.*, *J. Biol. Chem.* **283** (51), 35660-35667 (2008).

**Description:**

Human SMURF 1 (GenBank Accession No. NM\_020429), amino-acids 150-end, with N-terminal FLAG tag, MW=70 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

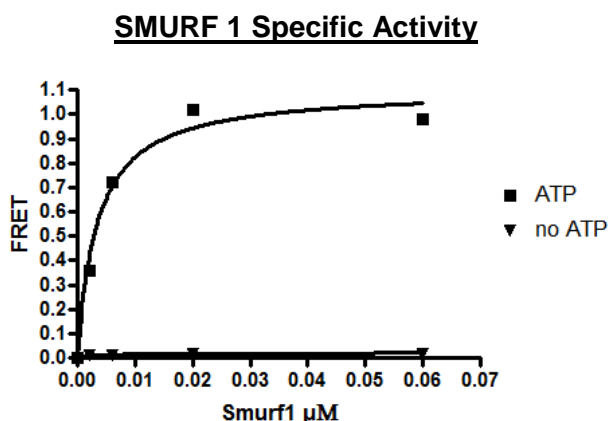
**Assay Conditions:**

10  $\mu\text{L}$  of reaction mixture contains 50 mM Tris, pH 7.4, 5 mM  $\text{MgCl}_2$ , 0.01% Tween, 1 mM DTT, 0.2  $\mu\text{M}$  Ub[Biotin], 50 nM Ube1, 1 mM ATP, 0.25  $\mu\text{M}$  of UBCH5b, and various amount of smurf1. The reaction was initiated with the addition of ATP, and incubated at  $30^{\circ}\text{C}$  for 3 hours. Tb labeled anti-Flag antibody donor and xl665 labeled acceptor was added and incubated at room temperature for hour.

**Application:**

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

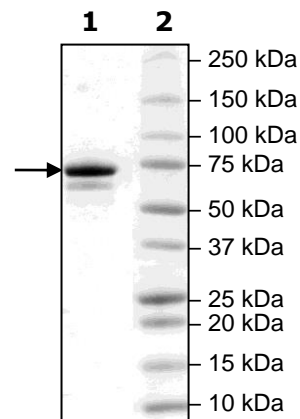
### Quality Assurance



**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2  $\mu\text{g}$  SMURF1  
**Lane 2:**  
Protein Marker

**MW:** 70 kDa  
**Purity:**  $\geq 85\%$



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)