

## Data Sheet

### UFL1, GST-tag

Human, recombinant, N-terminal GST-tag (CUL3)

**Catalog #:** 80411

**Lot #:** 150122      **Conc.:** 0.07 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.1% Tween-20, and 20% glycerol

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

#### References:

1. Lemaire, K. *et al.*, 2011. *PLOS One* **6(4)**: e18517.
2. Yang, P. *et al.*, 2014. *Immunobiol.* **219(6)**: 432-439.

**Description:** Human E3 UFM1-protein ligase 1 (UFL1), also known as LZAP-binding protein, GenBank Accession No. NM\_015323, a.a. 2-794 with N-terminal GST-tag, MW = 116 kDa, expressed in a baculovirus-infected Sf9 cell expression system.

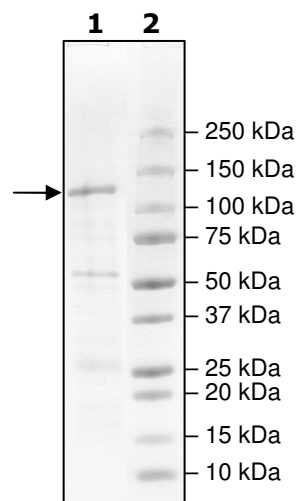
**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance

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

**Lane 1:**  
1.3  $\mu\text{g}$  UFL1  
**Lane 2:**  
Protein Marker

**MW:** 116 kDa  
**Purity:**  $\geq 55\%$



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)