

## Data Sheet

### UbchH7

Human, recombinant, C-terminal His tag

**Catalog #:** 80318

Lot #: 120103      Conc.: 0.89 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
 110 mM NaCl, 2.2 mM KCl, 200 mM  
 Imidazole, 3 mM DTT, and 20% glycerol.

**Stability:** >6 months at -80°C

**References:**

1. Katsanis, N., *et al.* 1998. *Genomics* **51**: 128-131.
2. Cook, W.J. *et al.*, 1997. *Biochemistry* **36** (7): 1621-1627.

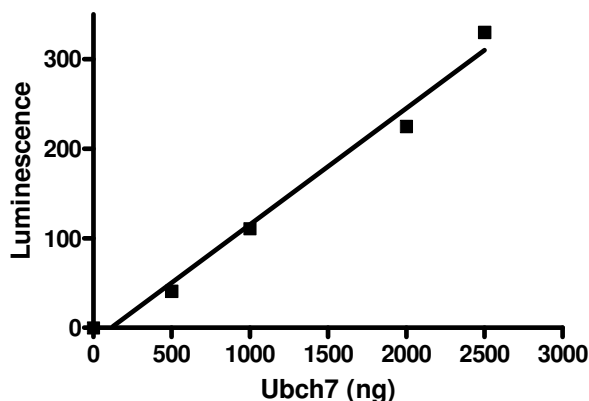
**Description:** Human UbchH7, also known as ubiquitin-conjugating enzyme E2L 3 (UBE2L3) (GenBank Accession No. NM\_003347), a.a. 1 – 154 (end) with C-terminal His tag, MW = 18.7 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Assay Conditions:** 50 µl reaction mix (4 mM ATP, 5 mM MgCl<sub>2</sub>, 1 mM DTT, UBE1 (0.5 µg), UbchH7, Biotin-Ubiquitin (0.5 µg) and Smurf2 (1.5 µg) in 50 mM Tris, pH 7.4) for 30 min at 37°C. Add to Streptavidin-coated plates for 30 min at 37°C. Incubate with anti-Smurf2, then HRP-labeled secondary antibody, and chemiluminescent detection.

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

## Quality Assurance

### Specific Activity

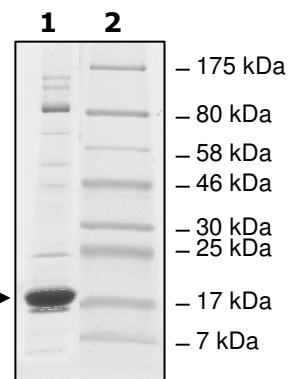


### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4.5 µg UbchH7

**Lane 2:**  
Protein Marker

**MW:** 18.7 kDa  
**Purity:** ≥71%



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)



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Sequence:**

**MAASRRLMKELEEIRKCGMKNFRNIQVDEANLLTWQGLIVPDNPPYDKGAFRIEINFP  
AEYPFKPPKITFKTKIYHPNIDEKGQVCLPVISAENWKPATKTDQVIQSLIALVNDPQPE  
HPLRADLAAEYSKDRKKFCKNAEEFTKKYGEKRPVDHHHHHH**

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)