

## Data Sheet

### UBA6 - UBE1L2

Human, recombinant, N-terminal Flag tag  
**Catalog #:** 80303  
**Lot #:** 100512                      **Conc:** 0.26mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
124 mM NaCl, 2.4 mM KCl, 90 µg/ml FLAG  
peptide, 10% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80°C

#### References:

1. Groettrup M. et al., Trends Biochem. Sci. 33 (5), 230-237 (2008).
2. Pelzer C. et al., J. Biol. Chem. 282 (32), 23010-23014 (2007).

#### Description:

Human UBA6 (UBE1L2) (GenBank Accession No. NM\_018227), amino-acids 2-end, with N-terminal Flag tag, MW=118 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

#### Application:

Useful in conjunction with E2 and E3 for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Assurance

### 10% SDS-PAGE Coomassie staining

