

## Data Sheet

### UBA6 - UBE1L2

Human, recombinant, N-terminal Flag tag  
**Catalog #:** 80303  
**Lot #:** 120523                      **Conc:** 0.11 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 0.005% Tween-  
20, 80 ng/ul Flag peptide, 20% Glycerol and  
1 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

