

Data Sheet

Ubch7 (UBE2L3)

Human, recombinant, N-terminal His-tag
Catalog #: 80317
Lot#91119 **Conc.** 0.89 mg/ml

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

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.

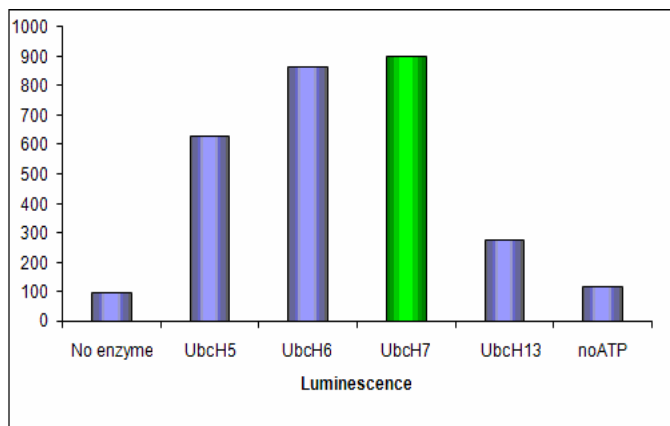
Description: Ubch7, also known as UBE2L3 (GenBank Accession No. NM_003347), amino acids 2-end, with N-terminal His-tag, MW = 18 kDa, expressed in an E. coli cell expression system.

Activity Assay: 50 µl reaction mix (4 mM ATP, 5 mM MgCl₂, 1 mM DTT, UBE1 (0.5 µg), Ubch7 or other E2, Biotin-Ubiquitin (0.5 µg) and Smurf2 (2 µg) in 50 mM Tris pH 7.4) for 30 min at 37°C, then add to streptavidin coated plates for 30 min at 37°C. Incubate with anti-Smurf2, then Europium-labeled secondary antibody, and detect by time resolved fluorescence (λ_{ex}=330 nm, λ_{em}=620 nm).

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

Quality Assurance

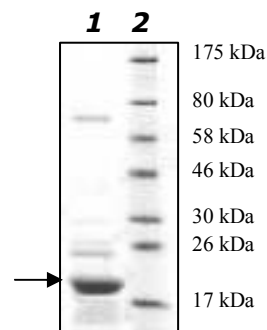
Ubch7 Activity



4-20 % SDS-PAGE Coomassie staining

Lane 1:
5 µg Ubch7
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 18 kDa
Purity: ~65%



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
 Please visit our website at: www.bpsbioscience.com