

## Data Sheet

### MCL1 His-Tag

Human, Recombinant, N-terminal His-Tag  
**Catalog #:** 40742  
**Lot #:** 180214      **Conc:** 0.8 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 0.04%  
Tween20, 3mM DTT, 20% glycerol,  
additional 200mM Imidazole

**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. Kozopas, K., *et al.*, *PNAS*. 1993; **90(8)**:  
3516-3520.
2. Leu, T., *et al.*, *New England Journal of  
Medicine*. 2013; **368**: 107-116.

#### Description:

Human MCL1, also known as Induced  
myeloid leukemia cell differentiation  
protein and BCL2L3, GenBank  
Accession No. NM\_021960, a.a. 2-327,  
with N-terminal His-tag, expressed in an  
E. coli expression system.

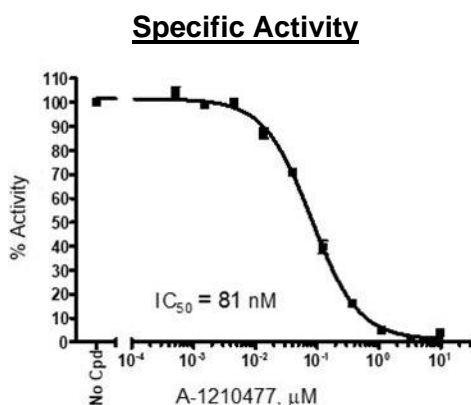
#### Assay Conditions:

Inhibition of MCL-1 by A-1210477, was  
measured using the MCL-1 TR-FRET  
Assay Kit, (BPS Bioscience #79506).

#### Applications:

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

## Quality Assurance



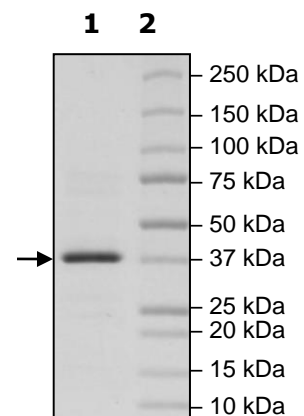
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
MCL1(2-327)

**Lane 2:**  
Protein Marker

**Purity:**  $\geq 90\%$

**MW:** 36 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.481.8694**, Fax  
**1.858.259.5734**

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)