

## Data Sheet

### IL-2 (C145A)

Human, Recombinant

**Catalog #:** 100159

**Lot #:** 180926

**Conc.:** 0.43 mg/ml

**Formulated in:** 10 mM Na phosphate  
pH7.4, 138 mM NaCl, 2.75 mM KCl.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid  
freeze/thaw cycles.

#### References:

1. Shevach EM. Trends Immunol. 2012 Dec,33(12):626-32.
2. Doms H, Abbas AK. Eur J Immunol. 2010 Jun,40(6):1538-40.

#### Description:

Human IL-2, also known as interleukin 2, (Genbank Accession No. NM\_000586), a.a. 21-153(end), with C145A mutation, expressed in an E.coli expression system. MW = 15 kDa.

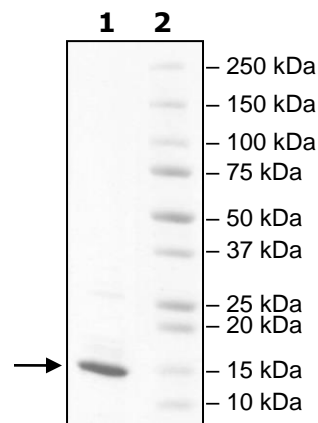
## Quality Assurance

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

**Lane 1:**  
2 $\mu\text{g}$  IL-2

**Lane 2:**  
Protein Marker

**MW:** 15 kDa  
**Purity:**  $\geq 90\%$



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)