

## Data Sheet

### PSMA, His-Avi-Tag

Human, Recombinant, N-terminal His-Avi-Tag  
**Catalog #:** 100463  
**Lot#:** 190820                      **Conc.:** 0.60 mg/ml

**Formulated in:** 8mM phosphate, pH 7.4,  
110 mM NaCl, 2.2mM KCl, and 20%  
glycerol.

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

#### References:

1. Hrkach, J. *et al.*, *Sci. Translational Med.* 2012; **4(128)**: 128-139.
2. Eiber, M. *et al.*, *J. Nucl. Med.* 2015; **56(5)**: 668-674.

**Description:** Human Prostate-specific membrane antigen (PSMA), also known as Folate Hydrolase 1, Glutamate Carboxypeptidase 2, FOLH, FGCP, and Cell Growth-Inhibiting Gene 27 Protein, GenBank Accession No. NM\_004476, a.a. 44-750(end) with an N-terminal His-Avi-Tag<sup>TM</sup>, expressed in a HEK293 cell expression system. MW = 84 kDa. This protein runs at a higher molecular weight due to glycosylation.

---

### Quality Assurance

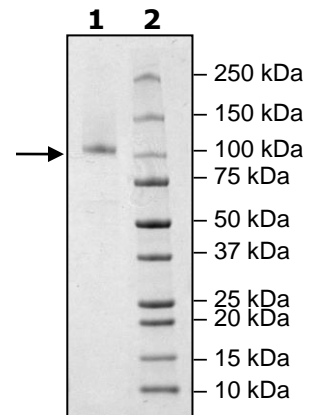
#### 4-20% SDS-PAGE Coomassie staining

Lane 1:  
4  $\mu\text{g}$  PSMA

Lane 2:  
Protein Marker

MW: 84 kDa+glycans

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 **1.858.202.1401** 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)