

## Data Sheet

### GPC3, Avi-His-Tagged

Human, Recombinant, C-terminal Avi-Tag™, C-terminal His-Tag

**Catalog #:** 100071

**Lot #:** 180508-1

**Conc.:** 0.82 mg/ml

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

**Stability:** At least 6 months at -80°C. Avoid freeze/thaw cycles. Protein may be diluted to ≥ 100 µg/ml in PBS + glycerol and stored at -80°C.

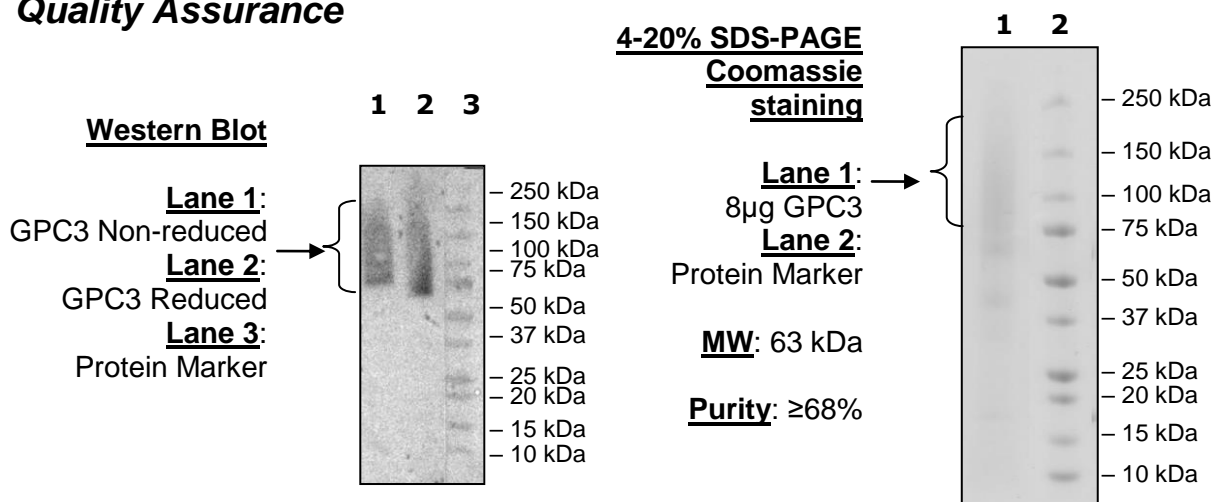
#### References:

1. Wykes, M., *et al.*, *Journal of Leukocyte Biology*, 2002 Oct; **72(4)**: 692-701.
2. Mukherjee, P., *et al.*, *Journal of Leukocyte Biology*. 2004 Jan; **77(1)**: 90-99.

#### Description:

Human GPC3, also known as Glypican 3, DGSX, and GTR2-2, with C-terminal Avi-Tag™ and C-terminal His-Tag. Genbank accession number NM\_004484, a.a. 25-559, expressed in a HEK293 cell expression system. MW = 63 kDa.

### Quality Assurance



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)