

## Data Sheet

### GPC3, Avi-His-Tag

Human, Recombinant, C-terminal Avi-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  $-80^{\circ}\text{C}$ . Avoid  
freeze/thaw cycles.

#### 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 Glypican 3, DGSX (GPC3), also known as GTR2-2, Genbank Accession No. NM\_004484, a.a. 25-559. with C-terminal Avi-His-tag™, expressed in a HEK293 cell expression system. MW = 63 kDa.

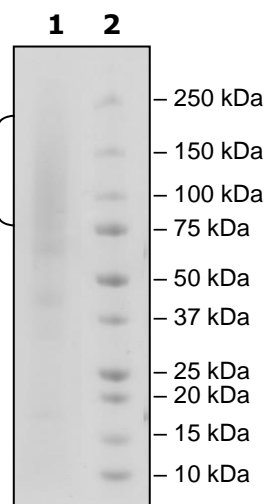
### Quality Assurance

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

**Lane 1:** →  
8µg GPC3  
**Lane 2:**  
Protein Marker

**MW:** 63 kDa

**Purity:** ≥68%



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)