

## Data Sheet

### **Mouse Sonic Hedgehog (Shh), His-tag**

Mouse, recombinant, C-terminal His-tag

**Catalog#:** 91012

**Lot#:** 170412A      **Conc.:** 0.57 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. Protein may be diluted to  $\geq 100\ \mu\text{g/ml}$  in PBS + glycerol and stored at  $-80^{\circ}\text{C}$ .

**Endotoxin Level:**  $<1\ \text{EU}/\mu\text{g}$

**References:**

1. Tran, P.V., *et al.*, *Nat. Genet.* 2008; **40**: 403-410.
2. Mukhopadhyay, S., *et al.*, *Cell* 2012; **152**(1-2): 210-223.

**Description:**

Mouse sonic hedgehog (mShh), also known as HHG-1. GenBank Accession No. NM\_009170, a.a. 25-198 with a C-terminal His-tag, expressed in an *E. coli* cell expression system, MW = 21 kDa.

**Application:**

Useful for studying protein binding and for screening small molecules and antibodies.

### **Quality Assurance**

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

**Lane 1:**

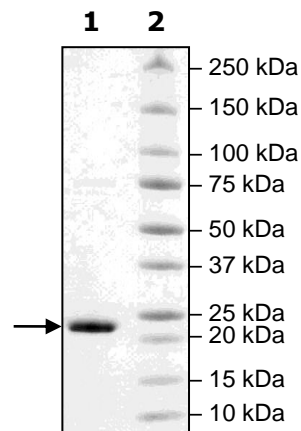
2  $\mu\text{g}$  Shh

**Lane 2:**

Protein Marker

**MW:** 21 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)