

## Data Sheet

### Luciferase, His-Tag (Firefly)

Firefly, Recombinant, C-terminal His-tag

**Catalog #:** 100576

**Lot #:** 200220-B

**Conc.:** 2.53 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 150 mM Imidazole and 20% glycerol.

**Description:** Firefly Luciferase, also known as Luciferin 4-monooxygenase, Uniprot No. Q27757, a.a. 1-545(full length) with C-terminal His-tag, expressed in an *E. coli* expression system. MW=61 kDa.

**Stability:** At least 6 months at -80°C. Avoid freeze/thaw cycles.

#### References:

1. DE Wet, J.R., *et al.*, *Molec. Cell. Bio.* 1987 Feb; **7(2)**:725-737.
2. Brasier, AR., *et al.*, *Euro. PMC.* 1989 Oct; **7(10)**:1116-1122.

---

### Quality Assurance

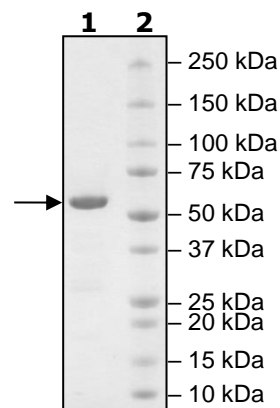
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2 µg Luciferase

**Lane 2:**  
Protein Marker

**MW:** 61 kDa

**Purity:** >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)