

## Data Sheet

### Sirtuin4

Human, recombinant, N-terminal GST tag  
**Catalog #:** 50015  
**Lot#:** 130812-Q2      **Conc:** 0.32 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
250 mM NaCl, and 20% glycerol.

**Stability:** >6 months at  $-70^{\circ}\text{C}$ . Avoid  
freeze/thaw cycles.

**References:**

1. Frye, R.A. (2000) *Biochem. Biophys. Res. Commun.* **273**, 793-798.

**Description:**

Human Sirtuin 4 (GenBank Accession No. NM\_012240), amino acids 25-314 with N-terminal GST tag, MW= 59 kDa, expressed in *E. coli* expression system.

**Application:**

Useful for screening inhibitors, and selectivity profiling.

## Quality Assurance

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

**Lane 1:**  
3  $\mu\text{g}$  SIRT4

**Lane 2:**  
Protein Marker

**MW:** 60 kDa  
**Purity:**  $\geq 85\%$

