

## Data Sheet

### HDAC3/NcoR2, His-Tag, GST-Tag

Human, recombinant, C-terminal His-tag,  
N-terminal GST-tag

**Catalog #:** 50003

**Lot#:** 190327

**Conc.:** 2.1 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 250 mM  
Imidazole and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid  
freeze/ thaw cycles.

#### References:

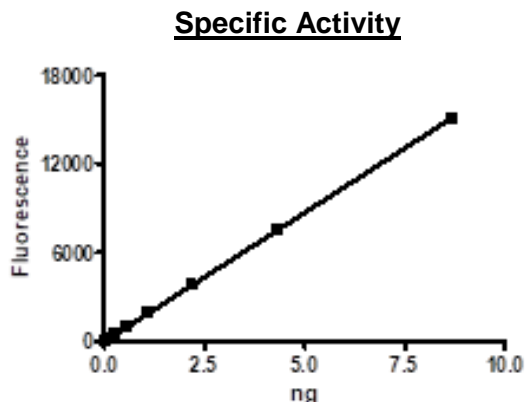
1. Gregoire, S. *et al. Mol. Cell. Biol.* **27** (4),  
1280-1295 (2007).
2. Escaffit, F. *et al., Mol. Cell. Biol.* **27** (2),  
554-567 (2007).

**Description:** Complex of human  
HDAC3, GenBank Accession No.  
NM\_003883, a.a. 1-428(end) with C-  
terminal His-tag, MW= 50 kDa, and  
human NCOR2, GenBank Accession  
No. NM\_006312, a.a. 395-489 with N-  
terminal GST tag, MW= 38 kDa, co-  
expressed in baculovirus expression  
system.

**Specific Activity:** 3000 pmol/min/ $\mu\text{g}$ .  
Assay conditions: 25 mM Tris/HCl, pH  
8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM  
MgCl<sub>2</sub>, and 0.1 mg/ml BSA, 20  $\mu\text{M}$  BPS  
HDAC substrate 3 (Catalog #50037),  
and HDAC3 (0.16 – 10ng). Incubation  
condition: 30 min at  $37^{\circ}\text{C}$  followed by  
developing for 15 min at room  
temperature.

**Application:** Useful for the study of  
enzyme kinetics, screening inhibitors,  
and selectivity profiling.

## Quality Assurance



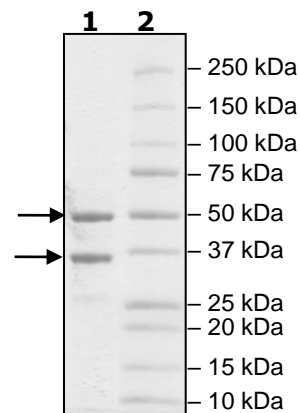
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4  $\mu\text{g}$  HDAC3/NcoR2

**Lane 2:**  
Protein Marker

**MW:** 50 kDa HDAC3  
38 kDa NcoR2

**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 **1.858.202.1401**, 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)