



11526 Sorrento Valley Rd SteA2  
San Diego Ca 92121  
Tel: 1.858.202.1401  
1.858.829.3082  
Fax: 1.858.259.5734  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### HDAC 3 / NcoR2

human, recombinant

Catalog #: 50003

Formulated in: 25 mM Tris-HCl, pH 8.0, 138 mM NaCl, 0.05% Tween-20, 20 mM glutathione and 10% glycerol.

Stability: >6 months at -80°C

References:

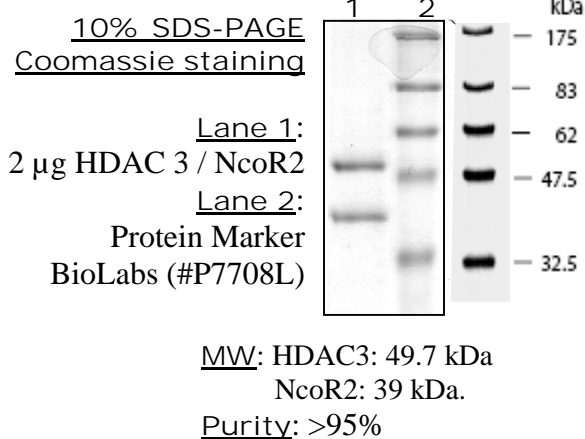
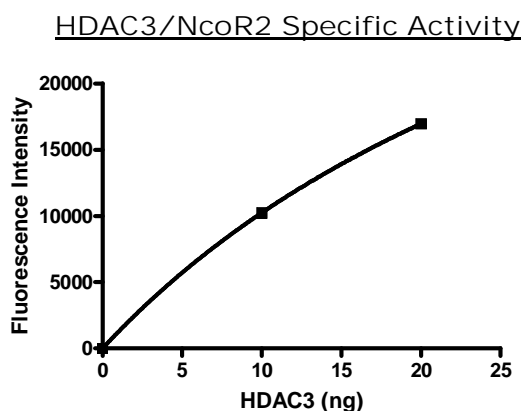
1. Gregoire, S. et al. Mol. Cell. Biol. 27 (4), 1280-1295 (2007)
2. Escaffit, F. et al., Mol. Cell. Biol. 27 (2), 554-567 (2007)

Description: Complex of human HDAC3 (GenBank Accession No. NM\_003883), full length with C-terminal His tag, MW= 49.7 kDa, and human NCOR2 (amino acid 395-489) (GenBank Accession No. NM\_006312), N-terminal GST tag, MW= 37.6 kDa, co-expressed in baculovirus expression system.

Specific Activity: 2500 pmol/min/μg. One U =1 pmol/min, Assay condition: 25 mM Tris/Cl, pH8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl<sub>2</sub>, and 0.1 mg/ml BSA, 30 μM BPS HDAC substrate (Catalog number 50032), and 1 ng/μl HDAC3/NcoR2. Incubation condition: 30 min at 37°C.

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

### Quality Assurance



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** or **1.858.829.3082**, Fax **1.858.259.5734**

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)



11526 Sorrento Valley Rd SteA2  
San Diego Ca 92121  
Tel: 1.858.202.1401  
1.858.829.3082  
Fax: 1.858.259.5734  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Assay Protocol**

**Material:** Assay buffer ( BPS catalog number 50031); Assay developer (BPS catalog number 50030); HDAC Substrate (BPS number 50032)

### **Step 1: adding all reaction mixture to a low binding black plate**

39 ul of HDAC assay buffer (BPS catalog number 50031)  
5 ul of 1 mg/ml BSA  
5 ul of 200 uM substrate (BPS catalog number: 50037)  
1 ul of HDAC3/NcoR2 (0.05 ug/ul)

Always add HDAC at the last.

Incubate at 37 °C for 30 min.

### **Setp2: stop the reaction**

add 50 ul of fluor de lys developer (2x) (BPS catalog number 50030) and incubate the plate at room temperature for 15 min

**Step 3:** read sample in a microtiter-plate reading fluorimeter capable of excitation at a wavelength in the range 350-380 nm and detection of emitted light in the range 440-460 nm.

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** or **1.858.829.3082**, Fax **1.858.259.5734**  
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)