



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Data Sheet**

### **HDAC Assay Developer**

**Catalog #: 50030**

**Lot#: 160922**

**DESCRIPTION:** HDAC assay developer is an essential reagent for signal generation in HDAC or sirtuin assays performed with any of BPS Bioscience's HDAC fluorimetric substrates (e.g. Cat. #50032). An appropriate fluorimetric substrate, comprising an acetylated lysine side chain, is first incubated with a sample containing HDAC activity (nuclear or cellular extract, purified enzyme, bead-bound immunocomplex, etc.). Deacetylation of the substrate sensitizes it so that, in a second step, treatment with the HDAC assay developer produces a fluorophore. Trichostatin A is included as an 'inhibitor stop' for class I and II HDACs.

**APPLICATIONS:** Study of HDAC or sirtuin kinetics, regulation and inhibitor sensitivity. Ideal for drug discovery and HTS applications and cell-based HDAC assays.

**SUPPLIED AS:** A 2x concentrate.

**QUANTITY:** 6 ml

**STORAGE:** -70°C. Store, long term, only as the 2x concentrate, at -70°C. Concentrate is stable for at least 3 freeze-thaw cycles, but care should be taken that both freezing and thawing steps are done quickly (e.g. liquid N<sub>2</sub> or dry ice/EtOH for freezing, rapid finger-rubbing for thawing). Frozen storage of smaller aliquots is recommended.

#### **REFERENCES:**

1. K. Ito *et al.* *Proc. Natl. Acad. Sci. USA* 2002 **99** 8921
2. B. Heltweg and M. Jung *Anal. Biochem.* 2002 **302** 175
3. S. Milutinovic *et al.* *J. Biol. Chem.* 2002 **277** 20974

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)