

## Data Sheet

### **pf-HDAC 1, His-FLAG-Tag**

*P. falciparum* HDAC1, recombinant,  
 C-terminal His-tag and FLAG-tag

**Catalog #:** 50063

**Lot #:** 100412      **Conc.:** 0.48 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
 124 mM NaCl, 2.4 mM KCl, 90 µg/ml FLAG  
 peptide, 10% glycerol.

**Stability:** >6 months at -80°C

**References:**

1. Chen *et al.*, *J. Med. Chem.* 2008 Jun 26;  
**51(12):**3437-48.
2. Joshi *et al.*, *Mol. Biochem. Parasitol.*  
 1999 Mar 15; **99(1):**11-9.

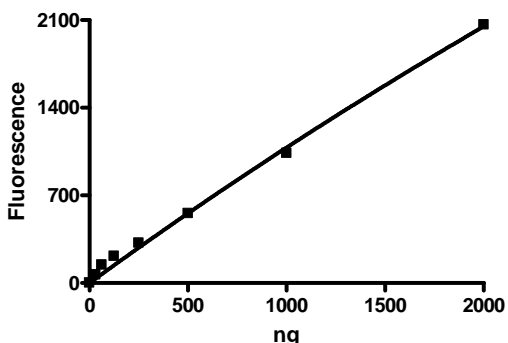
**Description:** Malarial (*Plasmodium falciparum*) HDAC, Genbank Accession No. AAD22407.1, full length (1-450) with C-terminal His-tag and FLAG-tag, MW=53 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Specific Activity:** 1.66 pmol/min/µg  
 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 µM BPS HDAC substrate 3 (Catalog #50037), and pf-HDAC1. HDAC reaction was done for 30 min at 37°C followed by treatment with 50 µl of HDAC developer (Cat. #50030) for 15 min at RT. Fluorescence intensity is measured at ex360/em460.

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

### Quality Assurance

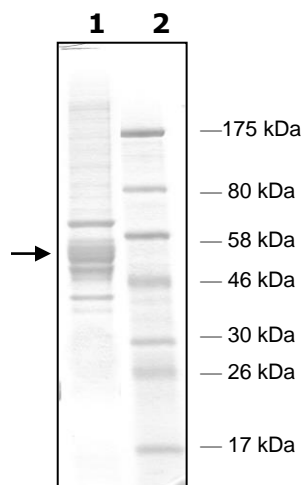
**Specific Activity**



**10% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
*pf*-HDAC 1  
**Lane 2:**  
 Protein Marker

**MW:** 53 kDa  
**Purity:** ≥58%



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)