

## Data Sheet

### HDAC11

Human, recombinant

**Catalog #:** 50021

**Lot #:** 141104

**Conc.:** 0.07 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 8.0, 500 mM NaCl, 50 mM KCl, 2 mM EDTA, 2 mM DTT, 0.5% Triton-X-100 and 10% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**Reference(s):**

1. Strausberg, R.L., *et al.*, *Proc. Natl. Acad. Sci. U.S.A.* **99** (26), 16899-16903 (2002).

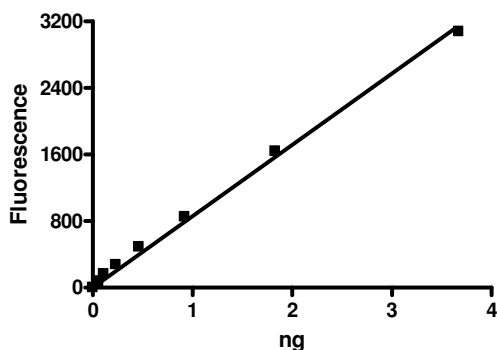
**Description:** Human HDAC11 (full length) (GenBank Accession No. NM\_024827) untagged, MW= 39.6 kDa, expressed in baculovirus expression system.

**Specific Activity:** 1500 pmol/min/ $\mu\text{g}$ .  
Assay conditions: 25 mM Tris/HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM  $\text{MgCl}_2$ , 0.05% Triton X-100, 0.1 mg/ml BSA, 20  $\mu\text{M}$  BPS HDAC substrate 1 (Catalog #50040), and HDAC11 (0.4 – 25 ng). 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

### Specific Activity

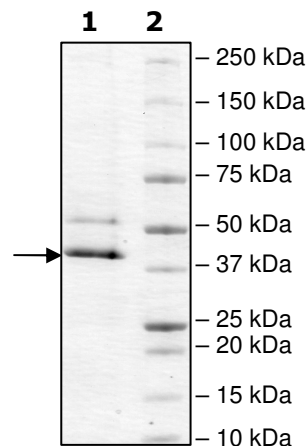


### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2  $\mu\text{g}$  HDAC11

**Lane 2:**  
Protein Marker

**MW:** 39.6 kDa  
**Purity:**  $\geq 70\%$



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)



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)

**Sequence:**

**GPLGSM**LHHTTQLYQHVPETRWPIVYSPRYNITFMGLEKLHPFDAGKWGKVINFLKEEKLLSDS  
MLVEAREASEEDLLVVHTRRYLNELKWSFAVATITEIPPVIFLPNFLVQRKVLRLRTQTGGTI  
MAGKLAVERGWAINVGGGFHHCSSDRGGGFCAYADITLAIKFLFERVEGISRATIIDLDAHQG  
NGHERDFMDDKRVIYIMDVYNRHIYPGDRFAKQAIRRKVELEWGTEDDEYLDKVERNIKKSLQ  
EHLPDVVVYNAGTDILEGDRLGGLSISPAGIVKRDELVFRMVRGRRVPILMVTSGGYQKRTAR  
IIADSILNLFGLGLIGPESPSVSAQNSDTPLLPPAVP

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)